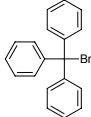
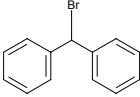
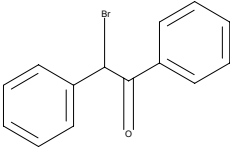
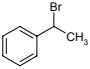
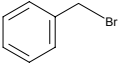
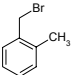
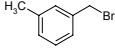
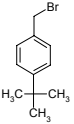
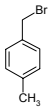
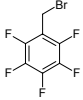
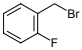
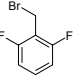
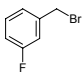
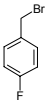
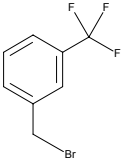
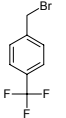
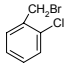
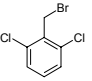
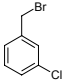
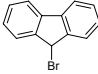


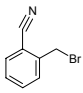
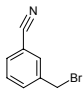
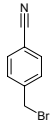
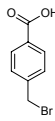
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD00000120	Brand ALDRICH	Product No. T83607	MDL MFCD00000134	Brand ALDRICH	Product No. B65403	MDL MFCD00000136	Brand ALDRICH	Product No. 258156	MDL MFCD00000139	Brand ALDRICH	Product No. 238104
formula $C_{19}H_{15}Br$	Molwgt 323.23	(C) 2008 SIGMA-ALDRICH CO.	formula $C_{13}H_{11}Br$	Molwgt 247.13	(C) 2008 SIGMA-ALDRICH CO.	formula $C_{14}H_{11}BrO$	Molwgt 275.14	(C) 2008 SIGMA-ALDRICH CO.	formula C_8H_9Br	Molwgt 185.06	(C) 2008 SIGMA-ALDRICH CO.
NAME Bromotriphenylmethane			NAME Bromodiphenylmethane			NAME Desyl bromide			NAME (1-Bromoethyl)benzene		
											

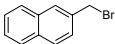
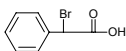
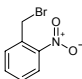
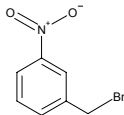
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD00000172	Brand ALDRICH	Product No. B17905	MDL MFCD00000174	Brand ALDRICH	Product No. B83401	MDL MFCD00000177	Brand ALDRICH	Product No. B83509	MDL MFCD00000180	Brand ALDRICH	Product No. 270768
formula C_7H_7Br	Molwgt 171.03	(C) 2008 SIGMA-ALDRICH CO.	formula C_8H_9Br	Molwgt 185.06	(C) 2008 SIGMA-ALDRICH CO.	formula C_8H_9Br	Molwgt 185.06	(C) 2008 SIGMA-ALDRICH CO.	formula $C_{11}H_{15}Br$	Molwgt 227.14	(C) 2008 SIGMA-ALDRICH CO.
NAME Benzyl bromide			NAME 2-Methylbenzyl bromide			NAME 3-Methylbenzyl bromide			NAME 4-tert-Butylbenzyl bromide		
											

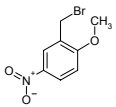
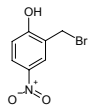
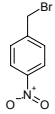
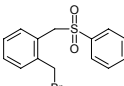
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD00000181	Brand DiscoveryCPR	Product No. CPR0181	MDL MFCD00000299	Brand ALDRICH	Product No. 101052	MDL MFCD00000324	Brand ALDRICH	Product No. 209511	MDL MFCD00000329	Brand ALDRICH	Product No. 264431
formula C_8H_9Br	Molwgt 185.06	(C) 2008 SIGMA-ALDRICH CO.	formula $C_7H_2BrF_5$	Molwgt 260.99	(C) 2008 SIGMA-ALDRICH CO.	formula C_7H_6BrF	Molwgt 189.02	(C) 2008 SIGMA-ALDRICH CO.	formula $C_7H_5BrF_2$	Molwgt 207.02	(C) 2008 SIGMA-ALDRICH CO.
NAME 4-METHYLBENZYL BROMIDE			NAME 2,3,4,5,6-Pentafluorobenzyl bromide			NAME 2-Fluorobenzyl bromide			NAME 2,6-Difluorobenzyl bromide		
											

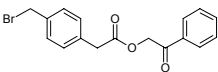
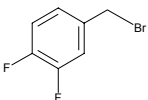
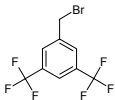
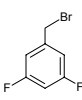
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD00000340	Brand ALDRICH	Product No. F7504	MDL MFCD00000359	Brand ALDRICH	Product No. 209538	MDL MFCD00000395	Brand ALDRICH	Product No. 290556	MDL MFCD00000403	Brand ALDRICH	Product No. 290564
formula C_7H_6BrF	Molwgt 189.02	(C) 2008 SIGMA-ALDRICH CO.	formula C_7H_6BrF	Molwgt 189.02	(C) 2008 SIGMA-ALDRICH CO.	formula $C_8H_6BrF_3$	Molwgt 239.03	(C) 2008 SIGMA-ALDRICH CO.	formula $C_8H_6BrF_3$	Molwgt 239.03	(C) 2008 SIGMA-ALDRICH CO.
NAME 3-Fluorobenzyl bromide			NAME 4-Fluorobenzyl bromide			NAME 3-(Trifluoromethyl)benzyl bromide			NAME 4-(Trifluoromethyl)benzyl bromide		
											

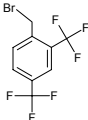
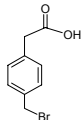
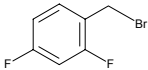
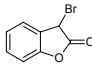
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD00000566	Brand ALDRICH	Product No. 252255	MDL MFCD00000577	Brand ALDRICH	Product No. 100307	MDL MFCD00000597	Brand ALDRICH	Product No. 136727	MDL MFCD00001133	Brand ALDRICH	Product No. B66604
formula C_7H_6BrCl	Molwgt 205.48	(C) 2008 SIGMA-ALDRICH CO.	formula $C_7H_5BrCl_2$	Molwgt 239.92	(C) 2008 SIGMA-ALDRICH CO.	formula C_7H_6BrCl	Molwgt 205.48	(C) 2008 SIGMA-ALDRICH CO.	formula $C_{13}H_9Br$	Molwgt 245.11	(C) 2008 SIGMA-ALDRICH CO.
NAME 2-Chlorobenzyl bromide			NAME 2,6-Dichlorobenzyl bromide - 97 %			NAME 3-Chlorobenzyl bromide			NAME 9-Bromofluorene		
											

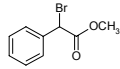
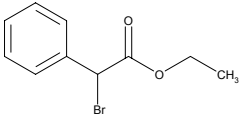
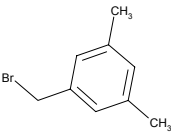
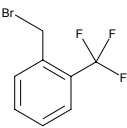
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD00001794	Brand ALDRICH	Product No. 144711	MDL MFCD00001809	Brand ALDRICH	Product No. 145610	MDL MFCD00001829	Brand ALDRICH	Product No. 144061	MDL MFCD00002567	Brand ALDRICH	Product No. 159549
formula C_8H_6BrN	Molwgt 196.04	(C) 2008 SIGMA-ALDRICH CO.	formula C_8H_6BrN	Molwgt 196.04	(C) 2008 SIGMA-ALDRICH CO.	formula C_8H_6BrN	Molwgt 196.04	(C) 2008 SIGMA-ALDRICH CO.	formula $C_8H_7BrO_2$	Molwgt 215.04	(C) 2008 SIGMA-ALDRICH CO.
NAME 2-(Bromomethyl)benzonitrile			NAME 3-(Bromomethyl)benzonitrile			NAME 4-(Bromomethyl)benzonitrile			NAME 4-(Bromomethyl)benzoic acid		
											

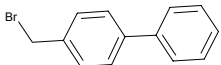
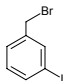
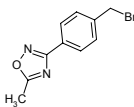
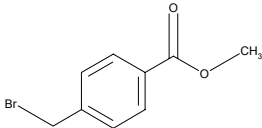
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD00004123	Brand ALDRICH	Product No. 143677	MDL MFCD00004206	Brand ALDRICH	Product No. B75859	MDL MFCD00007184	Brand DiscoveryCPR	Product No. CPR7184	MDL MFCD00007271	Brand ALDRICH	Product No. 222518
formula $C_{11}H_9Br$	Molwgt 221.09	(C) 2008 SIGMA-ALDRICH CO.	formula $C_8H_7BrO_2$	Molwgt 215.04	(C) 2008 SIGMA-ALDRICH CO.	formula $C_7H_6BrNO_2$	Molwgt 216.03	(C) 2008 SIGMA-ALDRICH CO.	formula $C_7H_6BrNO_2$	Molwgt 216.03	(C) 2008 SIGMA-ALDRICH CO.
NAME 2-(Bromomethyl)naphthalene			NAME alpha-Bromophenylacetic acid			NAME 2-NITROBENZYL BROMIDE			NAME 3-Nitrobenzyl bromide		
											

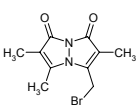
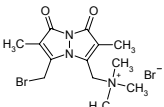
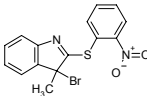
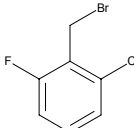
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD00007329	Brand ALDRICH	Product No. 163481	MDL MFCD00007340	Brand ALDRICH	Product No. 162906	MDL MFCD00007373	Brand ALDRICH	Product No. N13054	MDL MFCD00007552	Brand ALDRICH	Product No. 304123
formula $C_8H_8BrNO_3$	Molwgt 246.06	(C) 2008 SIGMA-ALDRICH CO.	formula $C_7H_6BrNO_3$	Molwgt 232.035	(C) 2008 SIGMA-ALDRICH CO.	formula $C_7H_6BrNO_2$	Molwgt 216.03	(C) 2008 SIGMA-ALDRICH CO.	formula $C_{14}H_{13}BrO_2S$	Molwgt 325.22	(C) 2008 SIGMA-ALDRICH CO.
NAME 2-Methoxy-5-nitrobenzyl bromide			NAME 2-Hydroxy-5-nitrobenzyl bromide			NAME 4-Nitrobenzyl bromide			NAME 1-Bromomethyl-2-((phenylsulfonyl)methyl)benzene		
											

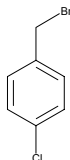
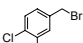
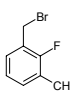
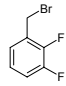
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD00009632	Brand ALDRICH	Product No. 317543	MDL MFCD00009889	Brand ALDRICH	Product No. 264458	MDL MFCD00009905	Brand ALDRICH	Product No. 263397	MDL MFCD00010304	Brand ALDRICH	Product No. 290319
formula $C_{17}H_{15}BrO_3$	Molwgt 347.2	(C) 2008 SIGMA-ALDRICH CO.	formula $C_7H_5BrF_2$	Molwgt 207.02	(C) 2008 SIGMA-ALDRICH CO.	formula $C_8H_5BrF_6$	Molwgt 307.03	(C) 2008 SIGMA-ALDRICH CO.	formula $C_7H_5BrF_2$	Molwgt 207.02	(C) 2008 SIGMA-ALDRICH CO.
NAME Phenacyl 4-(bromomethyl)phenylacetate			NAME 3,4-Difluorobenzyl bromide			NAME 3,5-Bis(trifluoromethyl)benzyl bromide			NAME 3,5-Difluorobenzyl bromide		
											

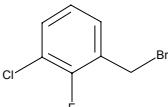
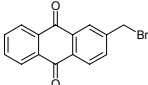
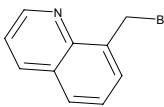
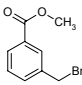
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD00010306	Brand ALDRICH	Product No. 290149	MDL MFCD00010632	Brand ALDRICH	Product No. 310417	MDL MFCD00011649	Brand ALDRICH	Product No. 264415	MDL MFCD00012320	Brand ALDRICH	Product No. 303453
formula $C_8H_5BrF_6$	Molwgt 307.03	(C) 2008 SIGMA-ALDRICH CO.	formula $C_9H_9BrO_2$	Molwgt 229.07	(C) 2008 SIGMA-ALDRICH CO.	formula $C_7H_5BrF_2$	Molwgt 207.02	(C) 2008 SIGMA-ALDRICH CO.	formula $C_8H_5BrO_2$	Molwgt 213.03	(C) 2008 SIGMA-ALDRICH CO.
NAME 2,4-Bis(trifluoromethyl)benzyl bromide			NAME 4-(Bromomethyl)phenylacetic acid, 97 %			NAME 2,4-Difluorobenzyl bromide			NAME 3-Bromo-2-coumaranone		
											

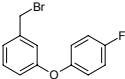
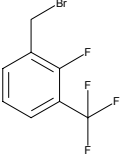
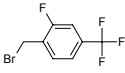
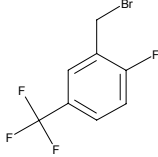
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD00013535	Brand ALDRICH	Product No. 365270	MDL MFCD00013536	Brand ALDRICH	Product No. 554065	MDL MFCD00013539	Brand DiscoveryCPR	Product No. CPS3539	MDL MFCD00013558	Brand ALDRICH	Product No. 254967
formula $C_9H_9BrO_2$	Molwgt 229.07	(C) 2008 SIGMA-ALDRICH CO.	formula $C_{10}H_{11}BrO_2$	Molwgt 243.1	(C) 2008 SIGMA-ALDRICH CO.	formula $C_9H_{11}Br$	Molwgt 199.09	(C) 2008 SIGMA-ALDRICH CO.	formula $C_8H_6BrF_3$	Molwgt 239.03	(C) 2008 SIGMA-ALDRICH CO.
NAME Methyl alpha-bromophenylacetate			NAME Ethyl alpha-bromophenylacetate			NAME 3,5-DIMETHYLBENZYL BROMIDE			NAME 2-(Trifluoromethyl)benzyl bromide		
											

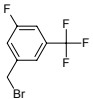
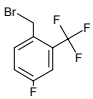
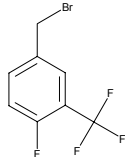
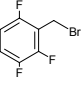
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD00017869	Brand DiscoveryCPR	Product No. CPS7869	MDL MFCD00019016	Brand ALDRICH	Product No. 427691	MDL MFCD00020808	Brand DiscoveryCPR	Product No. CPT0808	MDL MFCD00032453	Brand ALDRICH	Product No. 348155
formula $C_{13}H_{11}Br$	Molwgt 247.134	(C) 2008 SIGMA-ALDRICH CO.	formula C_7H_6BrI	Molwgt 296.93	(C) 2008 SIGMA-ALDRICH CO.	formula $C_{10}H_9BrN_2O$	Molwgt 253.098	(C) 2008 SIGMA-ALDRICH CO.	formula $C_9H_9BrO_2$	Molwgt 229.07	(C) 2008 SIGMA-ALDRICH CO.
NAME 4-BROMOMETHYLBIPHENYL			NAME 3-Iodobenzyl bromide			NAME 3-[4-(BROMOMETHYL)PHENYL]-5-METHYL-1,2,4-OXADIAZOLE			NAME Methyl 4-(bromomethyl)benzoate		
											

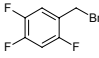
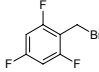
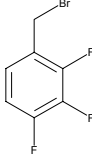
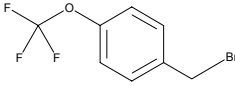
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD00036951	Brand SIAL	Product No. B4380	MDL MFCD00036953	Brand FLUKA	Product No. 71028	MDL MFCD00037978	Brand SIGMA	Product No. B4651	MDL MFCD00040126	Brand ALDRICH	Product No. 539090
formula $C_{10}H_{11}BrN_2O_2$	Molwgt 271.11	(C) 2008 SIGMA-ALDRICH CO.	formula $C_{13}H_{19}Br_2N_3O_2$	Molwgt 409.12	(C) 2008 SIGMA-ALDRICH CO.	formula $C_{15}H_{11}BrN_2O_2S$	Molwgt 363.23	(C) 2008 SIGMA-ALDRICH CO.	formula C_7H_5BrClF	Molwgt 223.47	(C) 2008 SIGMA-ALDRICH CO.
NAME Bromobimane			NAME Monobromo(trimethylammonio)bimane bromide			NAME 3-Bromo-3-methyl-2-(2-nitrophenylthio)-3H-indole			NAME 2-CHLORO-6-FLUOROBENZYL BROMIDE		
											

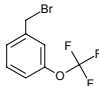
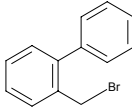
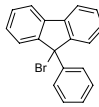
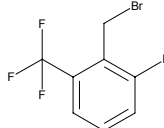
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD00040714	Brand DiscoveryCPR	Product No. CPV0714	MDL MFCD00041363	Brand ALDRICH	Product No. 681997	MDL MFCD00042470	Brand DiscoveryCPR	Product No. CPV2470	MDL MFCD00042488	Brand ALDRICH	Product No. 265314
formula C_7H_6BrCl	Molwgt 205.48	(C) 2008 SIGMA-ALDRICH CO.	formula $C_7H_5BrCl_2$	Molwgt 239.92	(C) 2008 SIGMA-ALDRICH CO.	formula C_8H_8BrF	Molwgt 203.053	(C) 2008 SIGMA-ALDRICH CO.	formula $C_7H_5BrF_2$	Molwgt 207.02	(C) 2008 SIGMA-ALDRICH CO.
NAME 4-CHLOROBENZYL BROMIDE			NAME 3,4-Dichlorobenzyl bromide			NAME 2-FLUORO-3-METHYLBENZYL BROMIDE, 95%			NAME 2,3-Difluorobenzyl bromide		
											

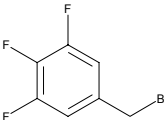
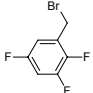
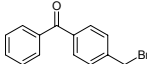
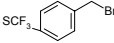
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD00042503	Brand DiscoveryCPR	Product No. CPV2503	MDL MFCD00042958	Brand ALDRICH	Product No. 444650	MDL MFCD00047621	Brand DiscoveryCPR	Product No. CPV7621	MDL MFCD00051437	Brand ALDRICH	Product No. 648116
formula C_7H_5BrClF	Molwgt 223.471	(C) 2008 SIGMA-ALDRICH CO.	formula $C_{15}H_9BrO_2$	Molwgt 301.13	(C) 2008 SIGMA-ALDRICH CO.	formula $C_{10}H_8BrN$	Molwgt 222.084	(C) 2008 SIGMA-ALDRICH CO.	formula $C_9H_9BrO_2$	Molwgt 229.07	(C) 2008 SIGMA-ALDRICH CO.
NAME 3-CHLORO-2-FLUOROBENZYL BROMIDE			NAME 2-Bromomethyl-antraquinone			NAME 8-(BROMOMETHYL)QUINOLINE			NAME Methyl 3-(bromomethyl)benzoate		
											

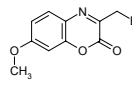
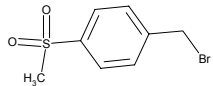
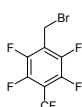
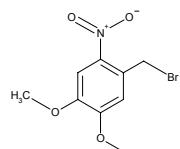
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD00059915	Brand DiscoveryCPR	Product No. CPW9915	MDL MFCD00061172	Brand DiscoveryCPR	Product No. CPX1172	MDL MFCD00061173	Brand DiscoveryCPR	Product No. CPX1173	MDL MFCD00061174	Brand DiscoveryCPR	Product No. CPX1174
formula C ₁₃ H ₁₀ BrFO	Molwgt 281.12	(C) 2008 SIGMA-ALDRICH CO.	formula C ₈ H ₅ BrF ₄	Molwgt 257.02	(C) 2008 SIGMA-ALDRICH CO.	formula C ₈ H ₅ BrF ₄	Molwgt 257.023	(C) 2008 SIGMA-ALDRICH CO.	formula C ₈ H ₅ BrF ₄	Molwgt 257.023	(C) 2008 SIGMA-ALDRICH CO.
NAME 3-(4-FLUOROPHENOXY)BENZYL BROMIDE			NAME 2-Fluoro-3-(trifluoromethyl)benzyl bromide			NAME 2-FLUORO-4-(TRIFLUOROMETHYL)BENZYL BROMIDE			NAME 2-FLUORO-5-(TRIFLUOROMETHYL)BENZYL BROMIDE		
											

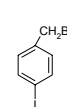
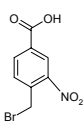
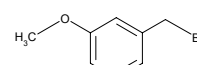
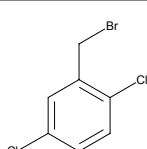
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD00061175	Brand DiscoveryCPR	Product No. CPX1175	MDL MFCD00061176	Brand ALDRICH	Product No. 436100	MDL MFCD00061177	Brand DiscoveryCPR	Product No. CPX1177	MDL MFCD00061208	Brand ALDRICH	Product No. 449407
formula C ₈ H ₅ BrF ₄	Molwgt 257.023	(C) 2008 SIGMA-ALDRICH CO.	formula C ₈ H ₅ BrF ₄	Molwgt 257.02	(C) 2008 SIGMA-ALDRICH CO.	formula C ₈ H ₅ BrF ₄	Molwgt 257.023	(C) 2008 SIGMA-ALDRICH CO.	formula C ₇ H ₄ BrF ₃	Molwgt 225.01	(C) 2008 SIGMA-ALDRICH CO.
NAME 3-FLUORO-5-(TRIFLUOROMETHYL)BENZYL BROMIDE			NAME 4-Fluoro-2-(trifluoromethyl)benzyl bromide			NAME 4-FLUORO-3-(TRIFLUOROMETHYL)BENZYL BROMIDE			NAME 2,3,6-Trifluorobenzyl bromide		
											

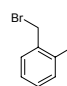
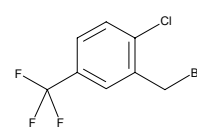
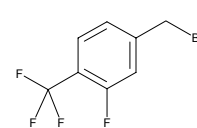
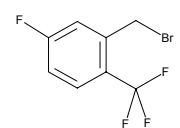
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD00061209	Brand DiscoveryCPR	Product No. CPX1209	MDL MFCD00061210	Brand DiscoveryCPR	Product No. CPX1210	MDL MFCD00061233	Brand ALDRICH	Product No. 554685	MDL MFCD00061238	Brand ALDRICH	Product No. 370622
formula C ₇ H ₄ BrF ₃	Molwgt 225.007	(C) 2008 SIGMA-ALDRICH CO.	formula C ₇ H ₄ BrF ₃	Molwgt 225.007	(C) 2008 SIGMA-ALDRICH CO.	formula C ₇ H ₄ BrF ₃	Molwgt 225.01	(C) 2008 SIGMA-ALDRICH CO.	formula C ₈ H ₆ BrF ₃ O	Molwgt 255.03	(C) 2008 SIGMA-ALDRICH CO.
NAME 2,4,5-TRIFLUOROBENZYL BROMIDE, 98%			NAME 2,4,6-TRIFLUOROBENZYL BROMIDE			NAME 2,3,4-Trifluorobenzyl bromide			NAME 4-(Trifluoromethoxy)benzyl bromide		
											

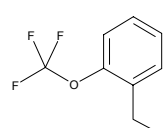
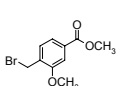
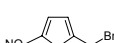
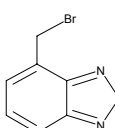
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD00061271	Brand ALDRICH	Product No. 408891	MDL MFCD00075454	Brand ALDRICH	Product No. 360082	MDL MFCD00075522	Brand ALDRICH	Product No. 368873	MDL MFCD00082477	Brand ALDRICH	Product No. 539627
formula C ₈ H ₆ BrF ₃ O	Molwgt 255.03	(C) 2008 SIGMA-ALDRICH CO.	formula C ₁₃ H ₁₁ Br	Molwgt 247.13	(C) 2008 SIGMA-ALDRICH CO.	formula C ₁₉ H ₁₃ Br	Molwgt 321.21	(C) 2008 SIGMA-ALDRICH CO.	formula C ₈ H ₅ BrF ₄	Molwgt 257.02	(C) 2008 SIGMA-ALDRICH CO.
NAME 3-(Trifluoromethoxy)benzyl bromide			NAME 2-Phenylbenzyl bromide			NAME 9-Bromo-9-phenylfluorene			NAME 2-Fluoro-6-(trifluoromethyl)benzyl bromide		
											

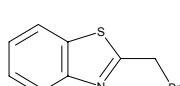
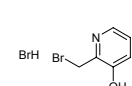
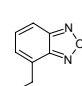
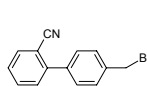
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD00083528	Brand ALDRICH	Product No. 559431	MDL MFCD00083543	Brand DiscoveryCPR	Product No. CPZ3543	MDL MFCD00092076	Brand ALDRICH	Product No. 449385	MDL MFCD00129189	Brand ALDRICH	Product No. 555444
formula C ₇ H ₄ BrF ₃	Molwgt 225.01	(C) 2008 SIGMA-ALDRICH CO.	formula C ₇ H ₄ BrF ₃	Molwgt 225.007	(C) 2008 SIGMA-ALDRICH CO.	formula C ₁₄ H ₁₁ BrO	Molwgt 275.14	(C) 2008 SIGMA-ALDRICH CO.	formula C ₈ H ₆ BrF ₃ S	Molwgt 271.1	(C) 2008 SIGMA-ALDRICH CO.
NAME 3,4,5-Trifluorobenzyl bromide			NAME 2,3,5-TRIFLUOROBENZYL BROMIDE			NAME 4-(Bromomethyl)benzophenone			NAME 4-(Trifluoromethylthio)benzyl bromide		
											

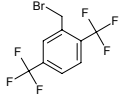
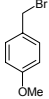
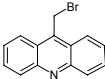
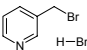
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD00143099	Brand FLUKA	Product No. 17631	MDL MFCD00185849	Brand DiscoveryCPR	Product No. CQJ5849	MDL MFCD00191855	Brand ALDRICH	Product No. 406406	MDL MFCD00192064	Brand ALDRICH	Product No. 392855
formula C ₁₀ H ₈ BrNO ₃	Molwgt 270.08	(C) 2008 SIGMA-ALDRICH CO.	formula C ₈ H ₉ BrO ₂ S	Molwgt 249.127	(C) 2008 SIGMA-ALDRICH CO.	formula C ₈ H ₂ BrF ₇	Molwgt 310.99	(C) 2008 SIGMA-ALDRICH CO.	formula C ₉ H ₁₀ BrNO ₄	Molwgt 276.08	(C) 2008 SIGMA-ALDRICH CO.
NAME 3-Bromomethyl-7-methoxy-1,4-benzoxazin-2-one			NAME 4-METHYLSULPHONYLBENZYL BROMIDE			NAME 2,3,5,6-Tetrafluoro-4-(trifluoromethyl)benzyl bromide			NAME 4,5-Dimethoxy-2-nitrobenzyl bromide		
											

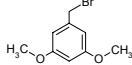
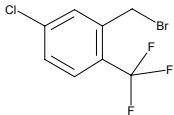
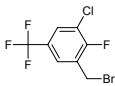
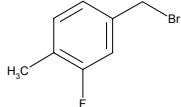
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD00209656	Brand ALDRICH	Product No. 515604	MDL MFCD00216575	Brand ALDRICH	Product No. 423564	MDL MFCD00216590	Brand ALDRICH	Product No. 429120	MDL MFCD00236034	Brand DiscoveryCPR	Product No. CQO6034
formula C ₇ H ₆ BrI	Molwgt 296.93	(C) 2008 SIGMA-ALDRICH CO.	formula C ₈ H ₆ BrNO ₄	Molwgt 260.04	(C) 2008 SIGMA-ALDRICH CO.	formula C ₈ H ₉ BrO	Molwgt 201.06	(C) 2008 SIGMA-ALDRICH CO.	formula C ₇ H ₅ BrCl ₂	Molwgt 239.92	(C) 2008 SIGMA-ALDRICH CO.
NAME 4-Iodobenzyl bromide			NAME 4-Bromomethyl-3-nitrobenzoic acid			NAME 3-Methoxybenzyl bromide			NAME 2,5-DICHLOROBENZYL BROMIDE		
											

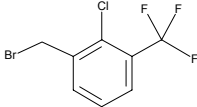
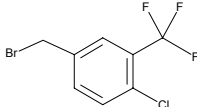
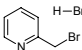
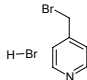
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD00236046	Brand ALDRICH	Product No. 634603	MDL MFCD00236216	Brand DiscoveryCPR	Product No. CQO6216	MDL MFCD00236289	Brand DiscoveryCPR	Product No. CQO6289	MDL MFCD00236290	Brand DiscoveryCPR	Product No. CQO6290
formula C ₇ H ₆ BrI	Molwgt 296.93	(C) 2008 SIGMA-ALDRICH CO.	formula C ₈ H ₅ BrClF ₃	Molwgt 273.478	(C) 2008 SIGMA-ALDRICH CO.	formula C ₈ H ₅ BrF ₄	Molwgt 257.023	(C) 2008 SIGMA-ALDRICH CO.	formula C ₈ H ₅ BrF ₄	Molwgt 257.023	(C) 2008 SIGMA-ALDRICH CO.
NAME 2-Iodobenzyl bromide			NAME 2-CHLORO-5-(TRIFLUOROMETHYL)BENZYL BROMIDE			NAME 3-FLUORO-4-(TRIFLUOROMETHYL)BENZYL BROMIDE			NAME 5-FLUORO-2-(TRIFLUOROMETHYL)BENZYL BROMIDE		
											

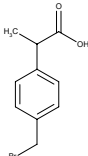
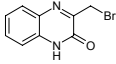
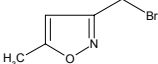
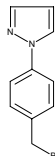
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD00236321	Brand DiscoveryCPR	Product No. CQO6321	MDL MFCD00270115	Brand ALDRICH	Product No. 681180	MDL MFCD00274235	Brand ALDRICH	Product No. 677329	MDL MFCD00457946	Brand DiscoveryCPR	Product No. CRK7946
formula C ₈ H ₆ BrF ₃ O	Molwgt 255.032	(C) 2008 SIGMA-ALDRICH CO.	formula C ₁₀ H ₁₁ BrO ₃	Molwgt 259.1	(C) 2008 SIGMA-ALDRICH CO.	formula C ₅ H ₄ BrNO ₃	Molwgt 205.99	(C) 2008 SIGMA-ALDRICH CO.	formula C ₇ H ₅ BrN ₂ S	Molwgt 229.1	(C) 2008 SIGMA-ALDRICH CO.
NAME 2-(TRIFLUOROMETHOXY)BENZYL BROMIDE			NAME Methyl 4-(bromomethyl)-3-methoxybenzoate			NAME 2-(Bromomethyl)-5-nitrofuran			NAME 4-(BROMOMETHYL)-2,1,3-BENZOTHIADIAZOLE		
											

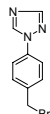
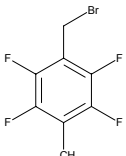
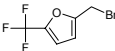
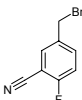
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD00461195	Brand DiscoveryCPR	Product No. CRL1195	MDL MFCD00466476	Brand DiscoveryCPR	Product No. CRL6476	MDL MFCD00515010	Brand DiscoveryCPR	Product No. CRQ5010	MDL MFCD00671503	Brand DiscoveryCPR	Product No. CSG1503
formula C ₈ H ₆ BrNS	Molwgt 228.112	(C) 2008 SIGMA-ALDRICH CO.	formula C ₆ H ₇ Br ₂ NO	Molwgt 268.935	(C) 2008 SIGMA-ALDRICH CO.	formula C ₇ H ₅ BrN ₂ O	Molwgt 213.033	(C) 2008 SIGMA-ALDRICH CO.	formula C ₁₄ H ₁₀ BrN	Molwgt 272.14	(C) 2008 SIGMA-ALDRICH CO.
NAME 2-(BROMOMETHYL)-1,3-BENZOTHIADIAZOLE			NAME 2-BROMOMETHYL-3-HYDROXYPYRIDINE HYDROBROMIDE			NAME 4-(BROMOMETHYL)-2,1,3-BENZOXADIAZOLE			NAME 4-BROMOMETHYL-2-CYANOBIIPHENYL		
											

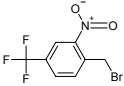
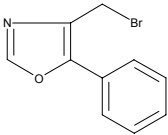
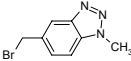
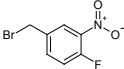
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD00799485	Brand ALDRICH	Product No. 465836	MDL MFCD00800380	Brand ALDRICH	Product No. 561282	MDL MFCD01321146	Brand DiscoveryCPR	Product No. CUT1146	MDL MFCD01321314	Brand ALDRICH	Product No. 475513
formula $C_9H_5BrF_6$	Molwgt 307.03	(C) 2008 SIGMA-ALDRICH CO.	formula C_8H_9BrO	Molwgt 201.06	(C) 2008 SIGMA-ALDRICH CO.	formula $C_{14}H_{10}BrN$	Molwgt 272.14	(C) 2008 SIGMA-ALDRICH CO.	formula $C_6H_7Br_2N$	Molwgt 252.93	(C) 2008 SIGMA-ALDRICH CO.
NAME 2,5-Bis(trifluoromethyl)benzyl bromide			NAME 4-Methoxybenzyl bromide			NAME 9-(Bromomethyl)acridine			NAME 3-(Bromomethyl)pyridine hydrobromide		
											

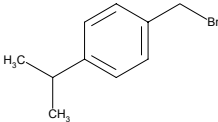
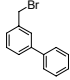
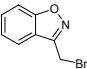
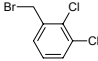
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD01321368	Brand ALDRICH	Product No. 480622	MDL MFCD01631356	Brand DiscoveryCPR	Product No. CVY1356	MDL MFCD01631513	Brand DiscoveryCPR	Product No. CVY1513	MDL MFCD01631534	Brand DiscoveryCPR	Product No. CVY1534
formula $C_9H_{11}BrO_2$	Molwgt 231.09	(C) 2008 SIGMA-ALDRICH CO.	formula $C_8H_5BrClF_3$	Molwgt 273.478	(C) 2008 SIGMA-ALDRICH CO.	formula $C_8H_4BrClF_4$	Molwgt 291.469	(C) 2008 SIGMA-ALDRICH CO.	formula C_8H_8BrF	Molwgt 203.053	(C) 2008 SIGMA-ALDRICH CO.
NAME 3,5-Dimethoxybenzyl bromide , 95 %			NAME 5-CHLORO-2-(TRIFLUOROMETHYL)BENZYL BROMIDE			NAME 3-CHLORO-2-FLUORO-5-(TRIFLUOROMETHYL)BENZYL BROMIDE			NAME 3-FLUORO-4-METHYLBENZYL BROMIDE		
											

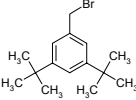
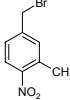
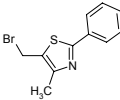
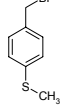
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD01631539	Brand DiscoveryCPR	Product No. CVY1539	MDL MFCD01631540	Brand DiscoveryCPR	Product No. CVY1540	MDL MFCD01863544	Brand ALDRICH	Product No. 491047	MDL MFCD01863545	Brand ALDRICH	Product No. 491748
formula $C_8H_5BrClF_3$	Molwgt 273.478	(C) 2008 SIGMA-ALDRICH CO.	formula $C_8H_5BrClF_3$	Molwgt 273.478	(C) 2008 SIGMA-ALDRICH CO.	formula $C_6H_7Br_2N$	Molwgt 252.93	(C) 2008 SIGMA-ALDRICH CO.	formula $C_6H_7Br_2N$	Molwgt 252.93	(C) 2008 SIGMA-ALDRICH CO.
NAME 2-CHLORO-3-(TRIFLUOROMETHYL)BENZYL BROMIDE			NAME 4-CHLORO-3-(TRIFLUOROMETHYL)BENZYL BROMIDE			NAME 2-(Bromomethyl)pyridine hydrobromide			NAME 4-(Bromomethyl)pyridine hydrobromide		
											

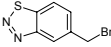
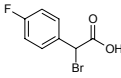
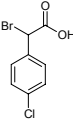
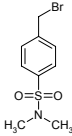
benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD02093445	Brand ALDRICH	Product No. 530360	MDL MFCD02660613	Brand ALDRICH	Product No. 534218	MDL MFCD02677720	Brand DiscoveryCPR	Product No. CZY7720	MDL MFCD02681940	Brand DiscoveryCPR	Product No. CZZ1940
formula $C_{10}H_{11}BrO_2$	Molwgt 243.1	(C) 2008 SIGMA-ALDRICH CO.	formula $C_9H_7BrN_2O$	Molwgt 239.07	(C) 2008 SIGMA-ALDRICH CO.	formula C_5H_5BrNO	Molwgt 176.012	(C) 2008 SIGMA-ALDRICH CO.	formula $C_{10}H_9BrN_2$	Molwgt 237.099	(C) 2008 SIGMA-ALDRICH CO.
NAME 2-[4-(Bromomethyl)phenyl]propanoic acid			NAME 3-(Bromomethyl)quinoxaline-2(1H)-one			NAME 3-(BROMOMETHYL)-5-METHYLISOXAZOLE			NAME 1-[4-(BROMOMETHYL)PHENYL]-1H-PYRAZOLE		
											

benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD02681943	Brand DiscoveryCPR	Product No. CZZ1943	MDL MFCD03001155	Brand ALDRICH	Product No. 556491	MDL MFCD03086219	Brand DiscoveryCPR	Product No. DBN6219	MDL MFCD03094341	Brand DiscoveryCPR	Product No. DBO4341
formula $C_9H_8BrN_3$	Molwgt 238.087	(C) 2008 SIGMA-ALDRICH CO.	formula $C_8H_5BrF_4$	Molwgt 257.02	(C) 2008 SIGMA-ALDRICH CO.	formula $C_6H_4BrF_3O$	Molwgt 228.995	(C) 2008 SIGMA-ALDRICH CO.	formula C_8H_5BrFN	Molwgt 214.036	(C) 2008 SIGMA-ALDRICH CO.
NAME 1-[4-(BROMOMETHYL)PHENYL]-1H-1,2,4-TRIAZOLE			NAME 4-Methyl-2,3,5,6-tetrafluorobenzyl bromide			NAME 2-(BROMOMETHYL)-5-(TRIFLUOROMETHYL)FURAN			NAME 3-CYANO-4-FLUOROBENZYL BROMIDE		
											

benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD03095413	Brand DiscoveryCPR	Product No. DBO5413	MDL MFCD03407316	Brand DiscoveryCPR	Product No. DCT7316	MDL MFCD03407338	Brand DiscoveryCPR	Product No. DCT7338	MDL MFCD03412241	Brand DiscoveryCPR	Product No. DCU2241
formula $C_8H_5BrF_3NO_2$	Molwgt 284.03	(C) 2008 SIGMA-ALDRICH CO.	formula $C_{10}H_8BrNO$	Molwgt 238.083	(C) 2008 SIGMA-ALDRICH CO.	formula $C_8H_8BrN_3$	Molwgt 226.076	(C) 2008 SIGMA-ALDRICH CO.	formula $C_7H_5BrFNO_2$	Molwgt 234.023	(C) 2008 SIGMA-ALDRICH CO.
NAME 2-NITRO-4-(TRIFLUOROMETHYL)BENZYL BROMIDE			NAME 4-(BROMOMETHYL)-5-PHENYL-1,3-OXAZOLE			NAME 5-(BROMOMETHYL)-1-METHYL-1H-1,2,3-BENZOTRIAZOLE			NAME 4-FLUORO-3-NITROBENZYL BROMIDE		
											

benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD03427264	Brand ALDRICH	Product No. 563285	MDL MFCD03452790	Brand ALDRICH	Product No. 538388	MDL MFCD03701148	Brand DiscoveryCPR	Product No. DDX1148	MDL MFCD03701317	Brand DiscoveryCPR	Product No. DDX1317
formula $C_{10}H_{13}Br$	Molwgt 213.11	(C) 2008 SIGMA-ALDRICH CO.	formula $C_{13}H_{11}Br$	Molwgt 247.13	(C) 2008 SIGMA-ALDRICH CO.	formula C_8H_8BrNO	Molwgt 212.045	(C) 2008 SIGMA-ALDRICH CO.	formula $C_7H_5BrCl_2$	Molwgt 239.926	(C) 2008 SIGMA-ALDRICH CO.
NAME 4-Isopropyl benzyl bromide			NAME 3-Phenylbenzyl bromide			NAME 3-(BROMOMETHYL)-1,2-BENZISOXAZOLE			NAME 2,3-DICHLOROBENZYL BROMIDE		
											

benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD03701623	Brand ALDRICH	Product No. 571474	MDL MFCD04039920	Brand ALDRICH	Product No. 572446	MDL MFCD04974048	Brand DiscoveryCPR	Product No. DIU4048	MDL MFCD05865196	Brand ALDRICH	Product No. 634816
formula $C_{15}H_{23}Br$	Molwgt 283.25	(C) 2008 SIGMA-ALDRICH CO.	formula $C_8H_8BrNO_2$	Molwgt 230.06	(C) 2008 SIGMA-ALDRICH CO.	formula $C_{11}H_{10}BNS$	Molwgt 268.177	(C) 2008 SIGMA-ALDRICH CO.	formula C_8H_9BrS	Molwgt 217.13	(C) 2008 SIGMA-ALDRICH CO.
NAME 3,5-Di-(tert-butyl)-benzyl bromide			NAME 3-Methyl-4-nitrobenzyl bromide			NAME 5-(BROMOMETHYL)-4-METHYL-2-PHENYL-1,3-THIAZOLE			NAME 4-(Methylthio)benzyl bromide		
											

benzyl bromides			benzyl bromides			benzyl bromides			benzyl bromides		
MDL MFCD06411539	Brand DiscoveryCPR	Product No. DOI1539	MDL MFCD08276759	Brand ALDRICH	Product No. 638668	MDL MFCD08276760	Brand ALDRICH	Product No. 638676	MDL MFCD08752694	Brand DiscoveryCPR	Product No. DXI2694
formula $C_7H_5BrN_2S$	Molwgt 229.1	(C) 2008 SIGMA-ALDRICH CO.	formula $C_8H_6BrFO_2$	Molwgt 233.03	(C) 2008 SIGMA-ALDRICH CO.	formula $C_8H_6BrClO_2$	Molwgt 249.49	(C) 2008 SIGMA-ALDRICH CO.	formula $C_9H_{12}BrNO_2S$	Molwgt 278.169	(C) 2008 SIGMA-ALDRICH CO.
NAME 5-(BROMOMETHYL)-1,2,3-BENZOTHIADIAZOLE			NAME alpha-Bromo-4-fluorophenylacetic acid			NAME alpha-Bromo-4-chlorophenylacetic acid			NAME 4-BROMOMETHYL-N,N-DIMETHYL-BENZENESULFONAMIDE		
											

benzyl bromides		
MDL MFCD09743698	Brand DiscoveryCPR	Product No. EBD3698
formula C_8H_8BrNO	Molwgt 214.061	(C) 2008 SIGMA-ALDRICH CO.
NAME 4-BROMOMETHYLBENZAMIDE		
