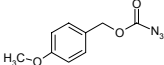
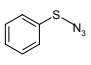
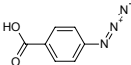
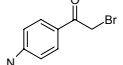
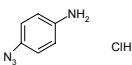
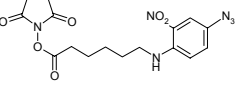
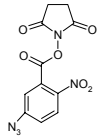
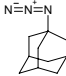
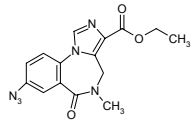
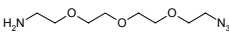
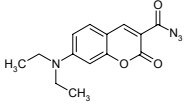
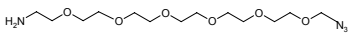


azides			azides			azides			azides		
MDL MFCD00001988	Brand ALDRICH	Product No. 152854	MDL MFCD00009700	Brand ALDRICH	Product No. 244546	MDL MFCD00013987	Brand DiscoveryCPR	Product No. CPS3987	MDL MFCD00042637	Brand FLUKA	Product No. 11550
formula C <sub>9</sub> H <sub>9</sub> N <sub>3</sub> O <sub>3</sub>	Molwgt 207.19	(C) 2008 SIGMA-ALDRICH CO.	formula C <sub>7</sub> H <sub>7</sub> N <sub>3</sub> S	Molwgt 165.22	(C) 2008 SIGMA-ALDRICH CO.	formula C <sub>7</sub> H <sub>5</sub> N <sub>3</sub> O <sub>2</sub>	Molwgt 163.135	(C) 2008 SIGMA-ALDRICH CO.	formula C <sub>8</sub> H <sub>6</sub> BrN <sub>3</sub> O	Molwgt 240.06	(C) 2008 SIGMA-ALDRICH CO.
NAME 4-Methoxybenzyloxycarbonyl azide			NAME Azidomethyl phenyl sulfide			NAME 4-AZIDOBENZOIC ACID, >98%			NAME 4-Azidophenacyl bromide		
											

azides			azides			azides			azides		
MDL MFCD00043274	Brand ALDRICH	Product No. 359556	MDL MFCD00049058	Brand SIGMA	Product No. A3407	MDL MFCD00054962	Brand SIGMA	Product No. A3282	MDL MFCD00074793	Brand SIAL	Product No. 276219
formula C <sub>6</sub> H <sub>7</sub> ClN <sub>4</sub>	Molwgt 170.6	(C) 2008 SIGMA-ALDRICH CO.	formula C <sub>16</sub> H <sub>18</sub> N <sub>6</sub> O <sub>6</sub>	Molwgt 390.35	(C) 2008 SIGMA-ALDRICH CO.	formula C <sub>11</sub> H <sub>7</sub> N <sub>5</sub> O <sub>6</sub>	Molwgt 305.2	(C) 2008 SIGMA-ALDRICH CO.	formula C <sub>10</sub> H <sub>15</sub> N <sub>3</sub>	Molwgt 177.25	(C) 2008 SIGMA-ALDRICH CO.
NAME 4-Azidoaniline hydrochloride			NAME 6-(4-Azido-2-nitrophenylamino)hexanoic acid N-hydroxysuccinimide ester			NAME 5-Azido-2-nitrobenzoic acid N-hydroxysuccinimide ester			NAME 1-Azidoadamantane		
											

azides			azides			azides			azides		
MDL MFCD00078598	Brand SIGMA	Product No. R109	MDL MFCD00269874	Brand FLUKA	Product No. 17758	MDL MFCD00467584	Brand FLUKA	Product No. 31755	MDL MFCD03453240	Brand FLUKA	Product No. 76172
formula C <sub>15</sub> H <sub>14</sub> N <sub>6</sub> O <sub>3</sub>	Molwgt 326.31	(C) 2008 SIGMA-ALDRICH CO.	formula C <sub>8</sub> H <sub>18</sub> N <sub>4</sub> O <sub>3</sub>	Molwgt 218.25	(C) 2008 SIGMA-ALDRICH CO.	formula C <sub>14</sub> H <sub>14</sub> N <sub>4</sub> O <sub>3</sub>	Molwgt 286.29	(C) 2008 SIGMA-ALDRICH CO.	formula C <sub>14</sub> H <sub>30</sub> N <sub>4</sub> O <sub>6</sub>	Molwgt 350.41	(C) 2008 SIGMA-ALDRICH CO.
NAME Ro 15-4513			NAME 11-Azido-3,6,9-trioxadecan-1-amine			NAME 7-(Diethylamino)coumarin-3-carbonyl azide			NAME O-(2-Aminoethyl)-O'-(2-azidoethyl)pentaethylene glycol		
											

azides		
MDL MFCD06798056	Brand ALDRICH	Product No. 573213
formula C <sub>14</sub> H <sub>19</sub> N <sub>3</sub> O <sub>3</sub>	Molwgt 277.32	(C) 2008 SIGMA-ALDRICH CO.
NAME (4S)-4-((1R)-2-Azido-1-(benzyloxy)ethyl)-2,2-dimethyl-1,3-dioxolane		
