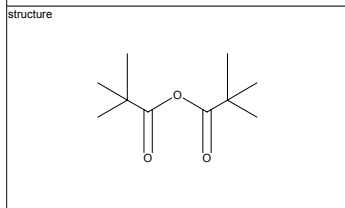


SIGMA -ALDRICH CO.

brand	Catalog number	formula	MW
Aldrich	143502	C ₁₀ H ₁₈ O ₃	186.25

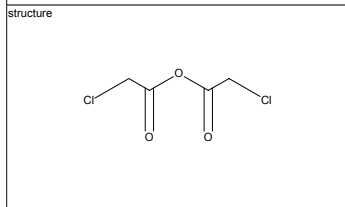
NAME
Pivalic anhydride, 99 %



SIGMA -ALDRICH CO.

brand	Catalog number	formula	MW
Aldrich	215163	C ₄ H ₄ Cl ₂ O ₃	170.98

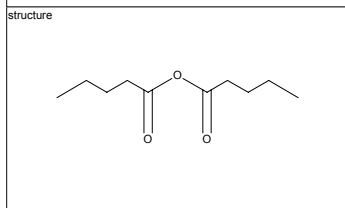
NAME
Chloroacetic anhydride, 97 %



SIGMA -ALDRICH CO.

brand	Catalog number	formula	MW
Aldrich	245933	C ₁₀ H ₁₈ O ₃	186.25

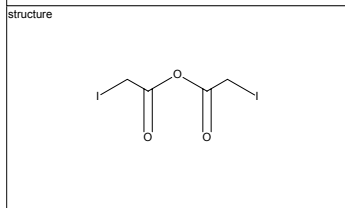
NAME
Valeric anhydride, 97 %



SIGMA -ALDRICH CO.

brand	Catalog number	formula	MW
Aldrich	284262	C ₄ H ₄ I ₂ O ₃	353.88

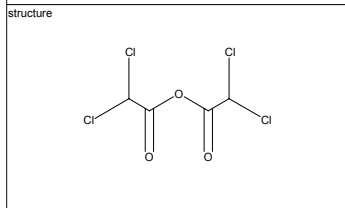
NAME
Iodoacetic anhydride



SIGMA -ALDRICH CO.

brand	Catalog number	formula	MW
Aldrich	331295	C ₄ H ₂ Cl ₄ O ₃	239.87

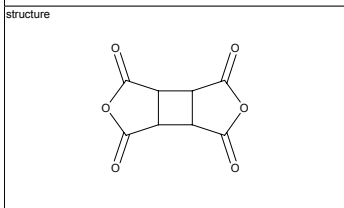
NAME
Dichloroacetic anhydride, gradetech. 85 %



SIGMA -ALDRICH CO.

brand	Catalog number	formula	MW
Aldrich	161330	C ₈ H ₄ O ₆	196.12

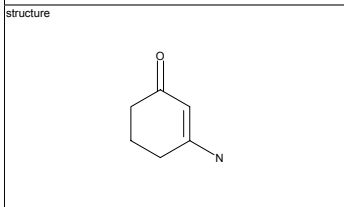
NAME
Cyclobutane-1,2,3,4-tetracarboxylic dianhydride



SIGMA -ALDRICH CO.

brand	Catalog number	formula	MW
Aldrich	239372	C ₆ H ₉ NO	111.14

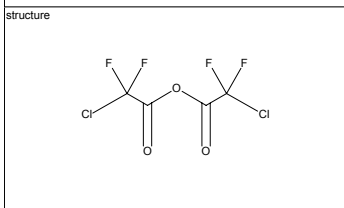
NAME
3-Amino-2-cyclohexen-1-one, 99 %



SIGMA -ALDRICH CO.

brand	Catalog number	formula	MW
Aldrich	252069	C ₄ Cl ₂ F ₄ O ₃	242.94

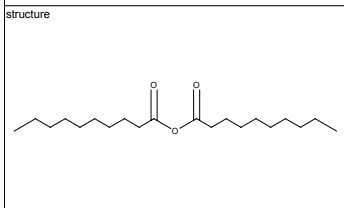
NAME
Chlorodifluoroacetic anhydride, 98 %



SIGMA -ALDRICH CO.

brand	Catalog number	formula	MW
Aldrich	284688	C ₂₀ H ₃₈ O ₃	326.52

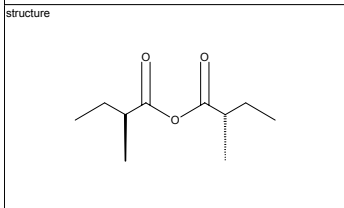
NAME
Decanoic anhydride, 98 %



SIGMA -ALDRICH CO.

brand	Catalog number	formula	MW
Aldrich	348562	C ₁₀ H ₁₈ O ₃	186.25

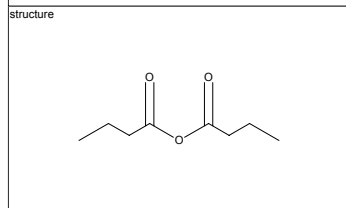
NAME
(S)-(+)-2-Methylbutyric anhydride, 94 %



SIGMA -ALDRICH CO.

brand	Catalog number	formula	MW
Aldrich	185736	C ₈ H ₁₄ O ₃	158.20

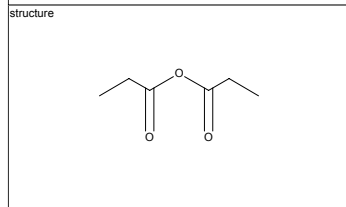
NAME
Butyric anhydride, 99 %



SIGMA -ALDRICH CO.

brand	Catalog number	formula	MW
Aldrich	240311	C ₆ H ₁₀ O ₃	130.14

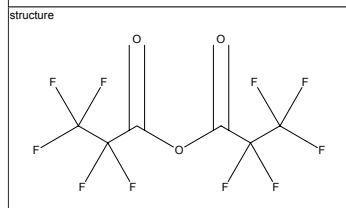
NAME
Propionic anhydride, 99+ %



SIGMA -ALDRICH CO.

brand	Catalog number	formula	MW
Aldrich	252387	C ₆ F ₁₀ O ₃	310.05

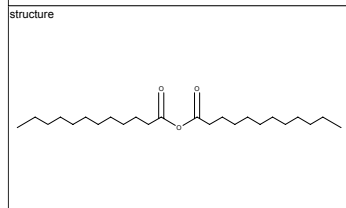
NAME
Pentafluoropropionic anhydride, 99 %



SIGMA -ALDRICH CO.

brand	Catalog number	formula	MW
Aldrich	286486	C ₂₄ H ₄₆ O ₃	382.63

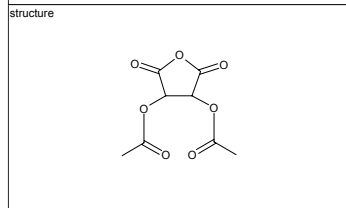
NAME
Dodecanoic anhydride, 98 %



SIGMA -ALDRICH CO.

brand	Catalog number	formula	MW
Aldrich	358924	C ₈ H ₈ O ₇	216.15

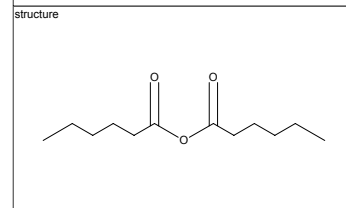
NAME
(+)-Di-O-acetyl-L-tartaric anhydride, 97 %



SIGMA -ALDRICH CO.

brand	Catalog number	formula	MW
Aldrich	194530	C ₁₂ H ₂₂ O ₃	214.31

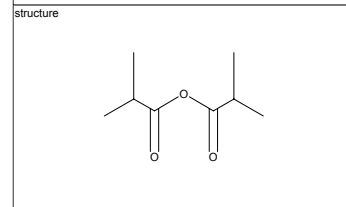
NAME
Hexanoic anhydride, 97 %



SIGMA -ALDRICH CO.

brand	Catalog number	formula	MW
Aldrich	245771	C ₈ H ₁₄ O ₃	158.20

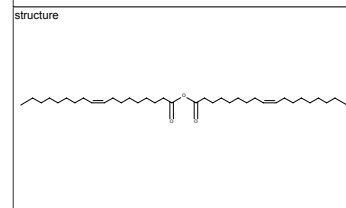
NAME
Isobutyric anhydride, 97 %



SIGMA -ALDRICH CO.

brand	Catalog number	formula	MW
Aldrich	284157	C ₃₆ H ₆₈ O ₃	546.93

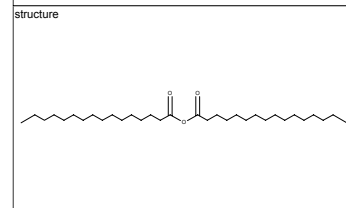
NAME
Oleic anhydride, 98 %



SIGMA -ALDRICH CO.

brand	Catalog number	formula	MW
Aldrich	286508	C ₃₂ H ₆₂ O ₃	494.85

NAME
Palmitic anhydride, 97 %



SIGMA -ALDRICH CO.

brand	Catalog number	formula	MW
Aldrich	367710	C ₄ Cl ₆ O ₃	308.76

NAME
Trichloroacetic anhydride, gradetech. 95 %

