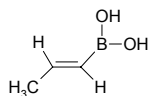


**Sigma - Aldrich
SAR Boronic Acids**

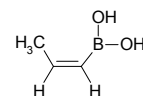
Brand	Catalog Number	MW
ALDRICH	165336	59.86

**Sigma - Aldrich
SAR Boronic Acids**

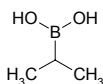
Brand	Catalog Number	MW
ALDRICH	576638	85.90

**Sigma - Aldrich
SAR Boronic Acids**

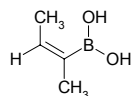
Brand	Catalog Number	MW
ALDRICH	572179	85.90

**Sigma - Aldrich
SAR Boronic Acids**

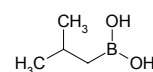
Brand	Catalog Number	MW
ALDRICH	648787	87.91

**Sigma - Aldrich
SAR Boronic Acids**

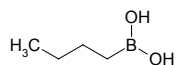
Brand	Catalog Number	MW
ALDRICH	638056	99.93

**Sigma - Aldrich
SAR Boronic Acids**

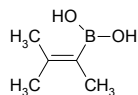
Brand	Catalog Number	MW
ALDRICH	346225	101.94

**Sigma - Aldrich
SAR Boronic Acids**

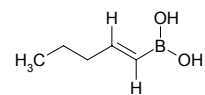
Brand	Catalog Number	MW
ALDRICH	163244	101.94

**Sigma - Aldrich
SAR Boronic Acids**

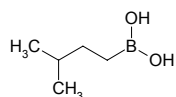
Brand	Catalog Number	MW
ALDRICH	639079	113.95

**Sigma - Aldrich
SAR Boronic Acids**

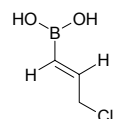
Brand	Catalog Number	MW
ALDRICH	578452	113.95

**Sigma - Aldrich
SAR Boronic Acids**

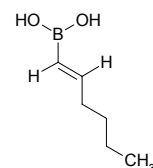
Brand	Catalog Number	MW
ALDRICH	655090	115.97

**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	556599	120.34

**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	521019	127.98





Sigma - Aldrich
SAR Boronic Acids

Brand	Catalog Number	MW
ALDRICH	521027	156.03

