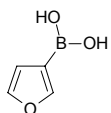




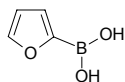
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	512168	111.89



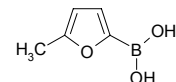
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	464910	111.89



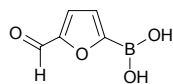
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	512265	125.92



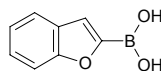
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	512346	139.90



**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	499943	161.95



**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	499951	212.01

