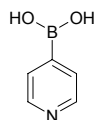




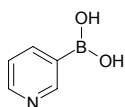
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	634492	122.92



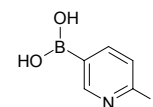
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	512125	122.92



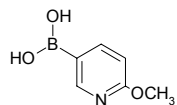
**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	639184	140.91



**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	637610	152.95



**Sigma - Aldrich
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	637386	157.37

