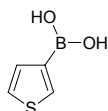




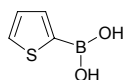
**Sigma - Aldrich  
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	436844	127.96



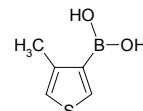
**Sigma - Aldrich  
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	436836	127.96



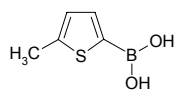
**Sigma - Aldrich  
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	542393	141.99



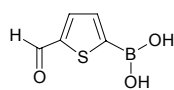
**Sigma - Aldrich  
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	512192	141.99



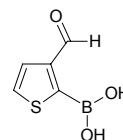
**Sigma - Aldrich  
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	514055	155.97



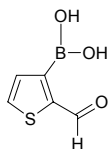
**Sigma - Aldrich  
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	499919	155.97



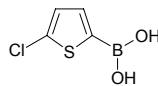
**Sigma - Aldrich  
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	499900	155.97



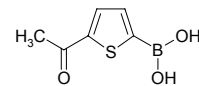
**Sigma - Aldrich  
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	499935	162.40



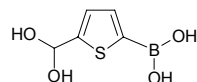
**Sigma - Aldrich  
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	499927	170.00



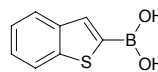
**Sigma - Aldrich  
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	470317	173.98



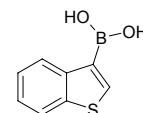
**Sigma - Aldrich  
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	499978	178.02



**Sigma - Aldrich  
SAR Boronic Acids**

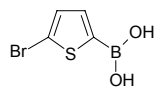
Brand	Catalog Number	MW
ALDRICH	512117	178.02





**Sigma - Aldrich**  
**SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	557684	206.85



**Sigma - Aldrich**  
**SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	499986	228.08

