



**Sigma - Aldrich  
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH		85.90



**Sigma - Aldrich  
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH		99.93



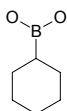
**Sigma - Aldrich  
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH		113.95



**Sigma - Aldrich  
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH		127.98



**Sigma - Aldrich  
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH		154.02

