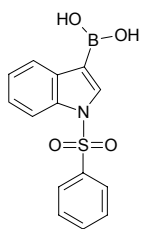




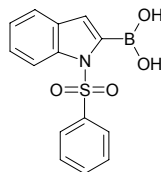
**Sigma - Aldrich  
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	563870	301.13



**Sigma - Aldrich  
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	563862	301.13



**Sigma - Aldrich  
SAR Boronic Acids**

Brand	Catalog Number	MW
ALDRICH	637394	339.98

