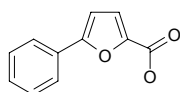




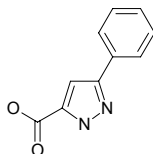
**Sigma - Aldrich**  
**SAR COOH**

Brand	Catalog Number	MW
ALDRICH	576301	188.18



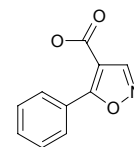
**Sigma - Aldrich**  
**SAR COOH**

Brand	Catalog Number	MW
ALDRICH	633739	188.19



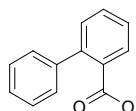
**Sigma - Aldrich**  
**SAR COOH**

Brand	Catalog Number	MW
ALDRICH	633682	189.17



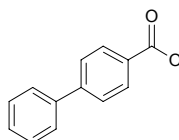
**Sigma - Aldrich**  
**SAR COOH**

Brand	Catalog Number	MW
ALDRICH	B34702	198.22



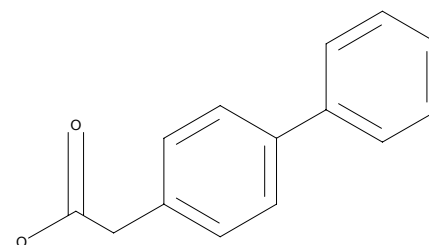
**Sigma - Aldrich**  
**SAR COOH**

Brand	Catalog Number	MW
ALDRICH	B34729	198.22



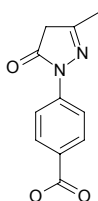
**Sigma - Aldrich**  
**SAR COOH**

Brand	Catalog Number	MW
ALDRICH	196487	212.25



**Sigma - Aldrich**  
**SAR COOH**

Brand	Catalog Number	MW
ALDRICH	197556	218.21



**Sigma - Aldrich**  
**SAR COOH**

Brand	Catalog Number	MW
ALDRICH	590827	228.25

