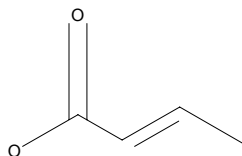




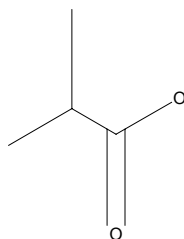
**Sigma - Aldrich**  
**SAR COOH**

Brand	Catalog Number	MW
ALDRICH	239569	86.09



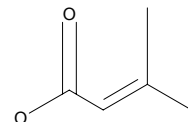
**Sigma - Aldrich**  
**SAR COOH**

Brand	Catalog Number	MW
ALDRICH	240168	88.11



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

Brand	Catalog Number	MW
ALDRICH	D138606	100.12



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

Brand	Catalog Number	MW
ALDRICH	277827	116.16

