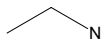




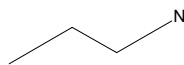
**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	395072	45.08



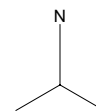
**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	240958	59.11



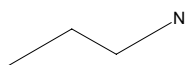
**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	109061	59.11



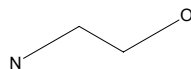
**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	109819	59.11



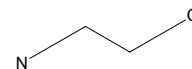
**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	411000	61.08



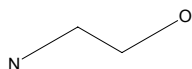
**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	110167	61.08



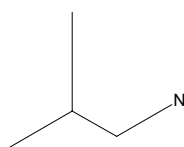
**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	398136	61.08



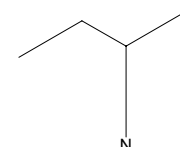
**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	114150	73.14



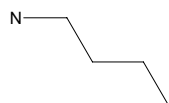
**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	B89000	73.14



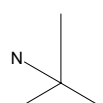
**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	471305	73.14



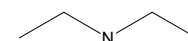
**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	391433	73.14



**Sigma - Aldrich  
SAR Amines**

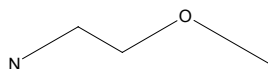
Brand	Catalog Number	MW
Aldrich	320277	73.14





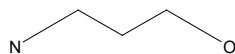
**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	241067	75.11



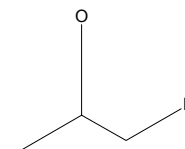
**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	239844	75.11



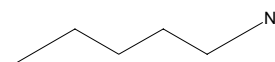
**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	110248	75.11



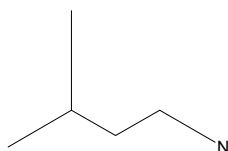
**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	171409	87.17



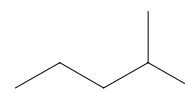
**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	126810	87.17



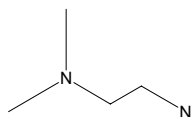
**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	171417	87.17



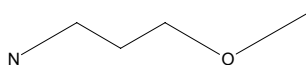
**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	D158003	88.15



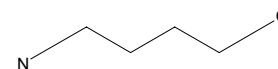
**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	M25007	89.14



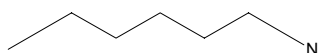
**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	178330	89.14



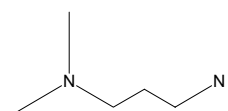
**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	219703	101.19



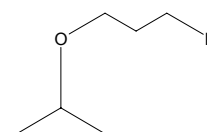
**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	240052	102.18



**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	123536	117.19





**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	O5802	129.25

