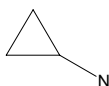




**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	125504	57.10



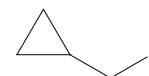
**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	394238	71.12



**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	359521	71.12



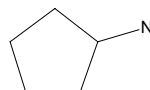
**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	225185	71.12



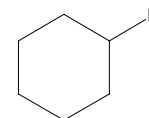
**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	C115002	85.15



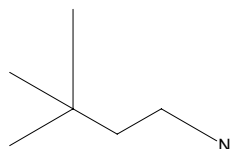
**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	240648	99.18



**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	183113	101.19



**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	159271	107.58

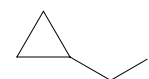


Cl



**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	A63805	107.58

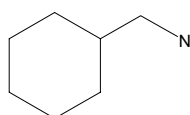


Cl



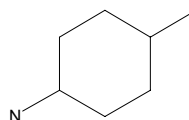
**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	101842	113.20



**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	177466	113.20



**Sigma - Aldrich  
SAR Amines**

Brand	Catalog Number	MW
Aldrich	255866	125.22

