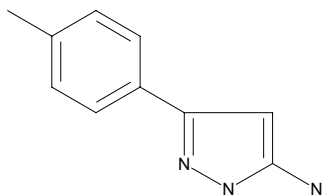




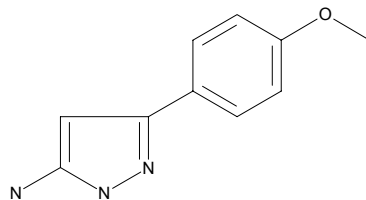
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	549207	173.22



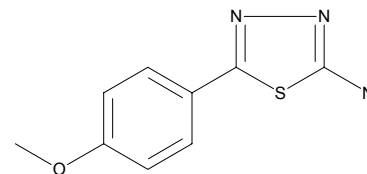
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	535370	189.22



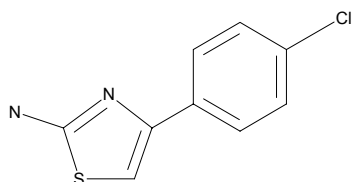
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	576026	207.26



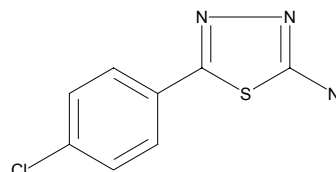
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	339369	210.69



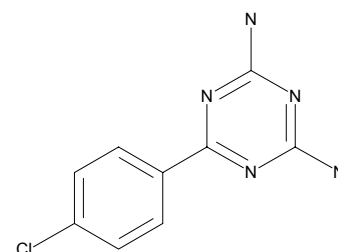
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	573744	211.67



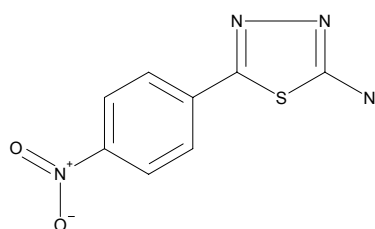
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	592188	221.65



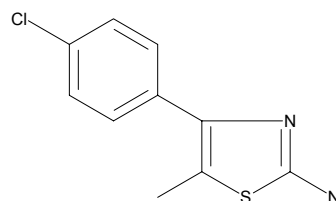
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	576018	222.23



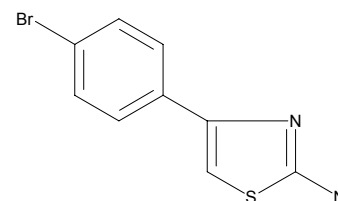
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
SALOR	L122033	224.71



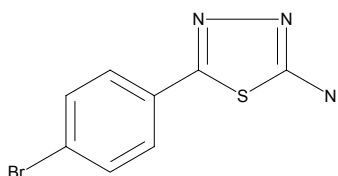
**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	534129	255.14



**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	576034	256.13



**Sigma - Aldrich
SAR Amines**

Brand	Catalog Number	MW
Aldrich	597244	266.10

