

Reagents for Surface Modification

Dojindo EU GmbH

Product Code	Product Name	Unit Size
Amino - Terminal		
A458-10	16-Amino-1-hexadecanethiol, hydrochloride	10 mg
A458-12		100 mg
A423-10	11-Amino-1-undecanethiol, hydrochloride	10 mg
A423-12		100 mg
A424-10	8-Amino-1-octanethiol, hydrochloride	10 mg
A424-12		100 mg
A425-10	6-Amino-1-hexanethiol, hydrochloride	10 mg
A425-12		100 mg
A483-10	Amino-EG ₆ -undecanethiol, hydrochloride	10 mg
A505-10	Amino-EG ₆ -hexadecanethiol, hydrochloride	10 mg
Amido - Terminal		
A508-10	5-Amido-pentanethiol	10 mg
A508-12		100 mg
A509-10	7-Amido-heptanethiol	10 mg
A509-12		100 mg
A510-10	10-Amido-decanethiol	10 mg
A510-12		100 mg
Fmoc-Amino-Terminal		
F287-10	N-Fmoc-Aminoundecanethiol	10 mg
F287-12		50 mg
F288-10	N-Fmoc-Aminooctanethiol	10 mg
F288-12		50 mg
F289-10	N-Fmoc-Aminoheptanethiol	10 mg
F289-12		50 mg
Carboxy - Terminal		
C488-10	Carboxylic acid-SAM Formation Reagent	1 ug x 3
C429-10	15-Carboxy-1-pentadecanethiol	10 mg
C385-10	10-Carboxy-1-decanethiol	10 mg
C385-12		100 mg
C386-10	7-Carboxy-1-heptanethiol	10 mg
C386-12		100 mg
C387-10	5-Carboxy-1-pentanethiol	10 mg
C387-12		100 mg
C445-10	Carboxy-EG ₆ -undecanethiol	10 mg
C445-12	Carboxy-EG ₆ -hexadecanethiol	100 mg
C463-10	Carboxy-EG ₆ -hexadecanethiol	10 mg
C404-10	10-Carboxydecyl disulfide	10 mg
C404-12		100 mg
C405-10	7-Carboxydecyl disulfide	10 mg
C406-10	5-Carboxydecyl disulfide	10 mg
D524-10	4,4'-Dithiodibutyric acid	500 mg

Product Code	Product Name	Unit Size
Hydroxy - Terminal		
H394-10	16-Hydroxy-1-hexadecanethiol	10 mg
H337-10		10 mg
H337-12	11-Hydroxy-1-undecanethiol	100 mg
H338-10		10 mg
H338-12	8-Hydroxy-1-octanethiol	100 mg
H354-10		10 mg
H354-12	Hydroxy-EG ₃ -undecanethiol	100 mg
H355-10		10 mg
H355-12	Hydroxy-EG ₆ -undecanethiol	100 mg
H395-10	Hydroxy-EG ₃ -hexadecanethiol	10 mg
H396-10	Hydroxy-EG ₆ -hexadecanethiol	10 mg
NHS ester-terminal		
D537-10		10 mg
D537-12	Dithiobis(succinimidyl undecanoate)	50 mg
D538-10		10 mg
D538-12	Dithiobis(succinimidyl octanoate)	50 mg
D539-10		10 mg
D539-12	Dithiobis(succinimidyl hexanoate)	50 mg
Ferrocenyl - Terminal		
F246-10		10 mg
F246-12	11-Ferrocenyl-1-undecanethiol	100 mg
F247-10		10 mg
F247-12	8-Ferrocenyl-1-octanethiol	100 mg
F269-10		10 mg
F269-12	6-Ferrocenyl-1-hexanethiol	100 mg
NTA - Terminal		
D550-10		10 mg
D550-12	Dithiobis(C ₂ -NTA)	50 mg
I279-10	Isothiocyanobenzyl-NTA	10 mg
A459-10	AB-NTA free acid	100 mg
M035-10	Maleimido-C ₃ -NTA	10 mg
EDTA- Terminal		
M030-10	Isothiocyanobenzyl-EDTA	10 mg
F279-10	FeBABE	1 mg
B437-10	BABE	10 mg
Sulfonate - Terminal		
S350-10	Sulfobetaine3-undecanethiol	10 mg
Biotin - Terminal		
B564-10	Biotin-SAM Formation Reagent	1 umol x 3
Carboxy - Terminal		
C488-10	Carboxylic acid-SAM Formation Reagent	1 umol x 3
Kit - Amine Coupling		
A515-10	Amine Coupling Kit	2 ml x 4