

SAMs 試薬 溶解性データ

品コード	製品 略名	Methyl alcohol (mol/l)				Chloroform (mol/l)				Ethyl alcohol (mol/l)				Dichloromethane(mol/l)				Tetrahydrofuran (mol/l)				Acetonitrile (mol/l)			
		10m	1m	100μ	10μ	10m	1m	100μ	10μ	10m	1m	100μ	10μ	10m	1m	100μ	10μ	10m	1m	100μ	10μ	10m	1m	100μ	10μ
A423	11-AUT,HCl													x	x			x	x	x		x	x	x	x
A424	8-AOT,HCl																					x			
A425	6-AHT,HCl					x								x				x				x	x	x	x
F287	11-FAUT	x								x												x			
F288	8-FAOT	x								x												x			
F289	6-FAHT																								
C385	10-CDT																								
C386	7-CHT																								
C387	5-CPT																								
C404	10-CDD					x								x				x				x			
C405	7-CHD																								
C406	5-CPD																								
D524	DDA													x								x			
D537	DSU																								
D538	DSO																								
D539	DSH																								
F246	11-FUT	x																							
F247	8-FOT																								
F269	6-FHT	x																							
H337	11-HUT																					x	x		
H338	8-HOT																								
H339	6-HHT																								
D550	Dithiobis(C2-NTA)					x	x	x	x					x	x	x	x	x	x	x	x	x	x	x	x

品コード	製品 略名	Ethyl acetate (mol/l)				Hexane (mol/l)				Dimethyl sulfoxide (mol/l)				Dimethyl formamide (mol/l)				Water (mol/l)			
		10m	1m	100μ	10μ	10m	1m	100μ	10μ	10m	1m	100μ	10μ	10m	1m	100μ	10μ	10m	1m	100μ	10μ
A423	11-AUT,HCl	x	x	x	x	x	x	x	x									x			
A424	8-AOT,HCl	x	x			x	x	x	x												
A425	6-AHT,HCl	x	x	x	x	x	x	x	x									x			
F287	11-FAUT					x	x	x	x									x	x	x	x
F288	8-FAOT					x	x	x	x									x	x	x	x
F289	6-FAHT					x	x	x	x									x	x	x	x
C385	10-CDT																	x	x	x	x
C386	7-CHT																	x			
C387	5-CPT																				
C404	10-CDD	x				x	x	x	x									x	x	x	x
C405	7-CHD					x	x	x	x									x	x	x	x
C406	5-CPD					x	x	x	x									x	x	x	x
D524	DDA					x	x	x	x									x	x	x	x
D537	DSU					x	x	x	x												
D538	DSO					x	x	x	x												
D539	DSH					x	x	x	x												
F246	11-FUT																	x	x	x	x
F247	8-FOT																	x	x	x	x
F269	6-FHT					x												x	x	x	x
H337	11-HUT					x												x	x	x	x
H338	8-HOT																	x			
H339	6-HHT																				
D550	Dithiobis(C2-NTA)	x	x	x	x	x	x	x	x	x				x							