thermoscientific



Choosing your SPE solution





Thermo Scientific[™] solid phase extraction (SPE) phases

Polymerics	Applications include
HyperSep [™] Retain PEP	Drugs and metabolites in biological matrices
Polystyrene divinylbenzene material surface modified with urea groups	i i i cita i o o o
HyperSep [™] Retain-CX Versatile polymeric material for retention of basic compounds	 Drugs of abuse from biological matrices HyperSep C18
HyperSep [™] Retain-AX Versatile polymeric material for retention of acidic compounds	 Acidic drugs of abuse from biologica matrices (THC and its metabolites)
HyperSep [™] Hypercarb Unique material for retention of	Retention and separation of highly polar species. Ideal for problem
highly polar compounds SOLA [™] and SOLAµ [™] HRP	 analytes in SPE applications Extraction of polar and non-polar analytes, such as vitamin D
Next-generation polystyrene divinylbenzene material surface functionalized with pyrrolidone	 biomarkers Drugs and metabolites in biological matrices Desalting of peptides in serum, plasma, or biological fluids
SOLA [™] and SOLAµ [™] SCX Next-generation polystyrene divinylbenzene material surface	 Enhanced retention of weak bases Drugs and metabolites in biological matrices, such as synthetic cathinones
functionalized with sulphonate groups	 Desalting of peptides in serum, plasma, or biological fluids
SOLA[™] and SOLAµ[™] SAX Next-generation polystyrene divinylbenzene material surface	 Enhanced retention of weak acids, such as 5-HIAA Drugs and metabolites in biological
functionalized with quaternary amine groups	 matrices Desalting of peptides in serum, plasma, or biological fluids
SOLA [™] and SOLAµ [™] WCX Next-generation polystyrene	 Enhanced retention of strong bases such as acetylcholinesterase inhibito Drugs and metabolites in biological
divinylbenzene material surface functionalized with carboxylic acid groups	 Desalting of peptides in serum, plasma, or biological fluids
SOLA [™] and SOLAµ [™] WAX Next-generation polystyrene	• Enhanced retention of strong acids, such as niflumic acid
divinylbenzene material surface functionalized with tertiary amine groups	 Drugs and metabolites in biological matrices Desalting of peptides in serum, plasma, or biological fluida.
Reversed-phase silica phases	plasma, or biological fluids Applications include
HyperSep [™] C18	Drugs and their metabolites in
Highly retentive alkyl-bonded silica phase for non-polar to moderately polar compounds	 biological matrices Trace organics in environmental wat samples Toxins in food samples
HyperSep [™] C8 Less retentive alternative to C18	 Drugs and their metabolites in biological matrices Trace organics in environmental wat
for non-polar to moderately polar compounds	samplesToxins in food samples
HyperSep [™] Phenyl Alternative selectivity for retention of basic compounds	Benzodiazepines in biologicalmatricesExtraction of aromatic compounds
Normal phase silica phases	Applications include
	AldehydesPesticides
HyperSep[™] Silica A polar sorbent primarily used to	CarotenoidsAflatoxins
retain analytes from non-polar	PhospholipidsAmines
matrices	HerbicidesFat soluble vitamins
HuporCop™ Flowie#	Fatty acids
HyperSep [™] Florisil Ideal for the isolation of polar	 Pesticides using AOAC and EPA methods, as well as
compounds from non-polar matrices	 Polychlorinated biphenyls (PCBs) in transformer oil
HyperSep[™] Cyano For retention of polar compounds from non-polar matrices	Retention of polar compounds from hexane and oil
HyperSep [™] Aminopropyl	Petroleum fractionation
A polar sorbent for both polar and anion exchange interactions	SaccharidesDrugs and drug metabolites
HyperSep[™] Diol For extraction of polar compounds	Normal phase extractionPurification of polar compounds
lon-exchange phases	Applications include
HyperSep [™] SAX (Strong anion exchanger) Strong anion exchange sorbent for extraction of weak acids	 Removal of acidic food pigments Removal of phenolic compounds Nucleic acids and surfactants
HyperSep [™] SCX	AntibioticsOrganic bases
(Strong cation exchanger) Strong cation exchange sorbent	Catecholamines
for extraction of charged basic compounds	DrugsAmino acidsHerbicides
HyperSep [™] Verify-CX Non-polar and anionic characteristics for improved analysis of basic drugs of abuse	 Basic drugs of abuse from biologica matrices

					non-polar solvent	moderately polar	polar solvent
Non-polar	Moderately polar	Polar	Anionic	Cationic	Non-polar	Moderately polar	Polar
Reversed-	Normal	Normal	Anion	Cation	Reversed-	Normal	Normal
phase	phase Reversed- phase	phase Reversed- phase	exchange	exchange	phase	phase Reversed- phase	phase
HyperSep Retain PEP	HyperSep Retain PEP	HyperSep Retain PEP	SOLA SAX	HyperSep Retain-CX	HyperSep Retain PEP	HyperSep Silica	HyperSep Hypercarb
SOLA HRP	SOLA HRP	HyperSep Hypercarb	SOLAµ SAX	SOLA SCX	SOLA HRP	HyperSep Florisil	HyperSep Cyano
SOLAµ HRP	SOLAµ HRP	HyperSep Cyano	SOLA WAX	SOLAµ SCX	SOLAµ HRP		HyperSep Aminopropy
HyperSep C18	HyperSep Silica	HyperSep Aminopropyl	SOLAµ WAX	SOLA WCX	HyperSep C18		HyperSep Diol
HyperSep C8	HyperSep Florisil	HyperSep Diol	HyperSep Retain-CX	SOLAµ WCX	HyperSep C8		
HyperSep Phenyl			HyperSep Retain-AX	HyperSep Verify-CX	HyperSep Phenyl		
			HyperSep Verify-AX	HyperSep SCX			
			HyperSep SAX				

HyperSep[™] Verify-AX Non-polar and cationic

Acidic drugs of abuse from biological matrices

characteristics for improved analysis of acidic drugs of abuse (THC and its metabolites)

Find out more at **thermofisher.com/speconsumables**

For Research Use Only. Not for use in diagnostic procedures. © 2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. This information is presented as an example of the capabilities of Thermo Fisher Scientific products. It is not intended to encourage use of these products in any manners that might infringe the intellectual property rights of others. Specifications, terms and pricing are subject to change. Not all products are available in all locations. Please consult your local sales representatives for details. **PO22067-EN 0620S**

ThermoFisher SCIENTIFIC