



revvity

For peace of mind, go NPE-Free.

Discover OptiScint™ NPE-Free Liquid Scintillation Solutions

Generate the high-quality data you need with our range of OptiScint NPE-free liquid scintillation cocktails and quenched standards.

Nonylphenol Ethoxylates (NPEs) are known to be toxic to humans, animals, and the environment. As a result, there are increasingly stringent regulations around products that contain NPEs.

While NPE-containing products, such as our Ultima Gold™ scintillation cocktails, can still be used for scientific R&D purposes under Article 3 of REACH regulations, some research labs want the option to use NPE-free products for greater peace of mind.



That's why Revvity has developed OptiScint cocktails and quenched standards, which do not contain NPEs either as a base or derivative component. This means they are compliant with regulations, such as the EEC directive 2003/53/EC and REACH, and you no longer need to worry about NPE-related safety concerns.

- No compromise on performance
- High flash point
- Safer handling
- Environmentally friendly

Product table for NPE-free optiscint cocktails

Part number	Product name	Description	Solvent	Flashpoint (°C)	Eff. without sample	Eff. with 10% water	Uptake capacity for water (mL)*
6013231 6013236	OptiScint MP OptiScint MP	NPE-Free Multipurpose Cocktail (1 L) NPE-Free Multipurpose Cocktail (5 L)	DIPN	>120	50	45	4
6013461 6013466	OptiScint LLT OptiScint LLT	NPE-Free Cocktail for Low Level (1 L) NPE-Free Cocktail for Low Level (5 L)	DIPN	116	46	40	10
6013281 6013286	OptiScint F OptiScint F	NPE-Free Cocktail for Filters (1 L) NPE-Free Cocktail for Filters (5 L)	DIPN	140	65	N/A	For organic samples and dried filter membranes only
6013791	OptiScint Flow OptiScint Flow	NPE-free Cocktail for Flow Scintillation Analyzers (1L) NPE-free Cocktail for Flow Scintillation Analyzers (5L)	DIPN	115	48	40	10

* This column shows the minimal uptake capacity for aqueous samples in 10 mL of cocktail.



| NPE-free OptiScint cocktails

For more information, visit www.revivity.com

