

Liquid scintillation counting with confidence.

Research and environmental labs around the world recognize that our liquid scintillation cocktails, tissue solubilizers, and specialty chemicals meet the highest standards of performance, convenience, and user and environmental safety – perfect for your most complex liquid scintillation counting challenges.

Combine our scintillation cocktails with trusted NEN® Radiochemicals, high-quality counting vials, market-leading radiometric detection instrumentation, and our portfolio of application notes for a complete single-source solution for your radiometric application needs.

#### Safer scintillation cocktails

#### Ultima Gold family

Ultima Gold™ is a safer liquid scintillation cocktail for a wide range of aqueous and nonaqueous samples. This multipurpose LSC cocktail has a high counting efficiency and provides superior detection for samples that exhibit severe quench in conventional cocktails



Product	Product no.	Size
	6013326	1 x 5 L
<b>Ultima Gold</b> - Safe scintillation cocktail for wide range of aqueous and nonaqueous samples	6013327	4 x 2.5 L
Horiaqueous samples	6013329	2 x 5 L
Ultima Gold F - For counting dry filters and nonaqueous (organic) samples	6013179	2 x 5 L
Illtima Cald VD. Dravidas vary bigh savarala load savarity.	6013117	4 x 2.5 L
Ultima Gold XR - Provides very high sample-load capacity	6013119	2 x 5 L
Ultima Gold MV - For rapid uptake samples, counting wet glass fiber filters	6013151	1 x 1 L
from cell harvesters	6013159	2 x 5 L
<b>Ultima Gold LLT</b> - Low levels of <sup>3</sup> H in wide-range water up to 54%, counting efficiencies of 30%		1 x 5 L
		4 x 2.5 L
<b>Ultima Gold uLLT -</b> for ultra low level LSC applications. The ultra low background, high efficiency and good sample uptake is optimal for	6013681	1 x 1 L
environmental monitoring applications.	6013687	4 x 2.5 L
Ultima Gold AB - Alpha/beta discrimination with slow pulse-decay	6013301	2 x 5 L
characteristics	6013309	2 x 5 L
OptiPhase HiSafe 3 - High counting efficiency for nonaqueous samples	1200-437	1 x 5 L
<b>Emulsifier-Safe</b> - LAB-based for aqueous and organic samples up to 15% sample load	6013389	2 x 5 L

#### Classical cocktails

Our classical liquid scintillation cocktail formulations are optimized for the highest counting efficiency and maximum sample-holding capacity. Prepared cocktails are easy to use, save prep time, and minimize errors. And our carefully controlled blending and quality-assurance procedures provide high performance, batch homogeneity, and lot-to-lot uniformity.

Product	Product no.	Size
<b>BetaPlate Scint</b> - For samples harvested or spotted on dry filter membranes; ideal for organic samples	1205-440	1 x 5 L
High-Efficiency Mineral Oil Scintillator - For detecting radon in water and soil samples	6NE9571	1 x 1 L
Insta-Gel Plus - Ideal for counting large	6013391	1 x 1 L
volumes of water with suspended solids	6013399	2 x 5 L
<b>Pico-Fluor Plus</b> - ideal for counting aqueous samples and buffer solutions	6013691	1 x 1 L
and is compatible with alkaline tissue solubilizers and alkaline samples.	6013699	2 x 5 L
Insta-Fluor Plus	6013167	4 x 2.5 L
Hionic-Fluor - ideal for counting	6013311	1 x 1 L
alkaline and concentrated salt samples.	6013319	2 x 5 L

#### Cocktails for microplate formats

Our MicroScint™ cocktails mix easily with samples for rapid uptake of aqueous and nonaqueous samples in microplates. These safer cocktails provide you with a choice of optimal characteristics, including excellent counting efficiency, high sample capacity, quench resistance, and polystyrene compatibility.

Product	Product no.	Size
<b>MicroScint-20</b> - Accepts diluted aqueous samples up to 20% loading, damp filters	6013621	1 x 5 L
<b>MicroScint-O</b> - For counting nonpolar organic samples for dry filters	6013611	1 x 5 L
<b>MicroScint-PS</b> - Formulated for polystyrene; high counting stability, up to 20% loading	6013631	2 x 5 L
MetiLex Solid Scintillator - For Microbeta with filtermat harvested	1450-441	100/pack
or dot-blotted samples	1450-442	75/pack

#### Cocktails for filters

Filter Count will dissolve these wet or dry filters, reducing sample preparation.

Product	Product no.	Size
Filter-Count™ - used for counting wet	6013141	1 x 1 L
and dry membrane filters.	6013149	2 x 5 L

#### Flow-detection cocktails

The high loading capacity of Ultima-Flo™ Flo-Scint® means less cocktail required, longer residence times, and higher sensitivity.

Product	Product no.	Size
Ultima-Flo AP - Flow counting;	6013591	1 x 1L
accepts gradients up to 2.0 M ammonium phosphate	6013599	2 x 5 L
Flo-Scint II	6013529	2 x 5L
Flo-Scint A	6013569	2 x 5L

# Sample oxidizer cocktails

These pseudocumene-based oxidizer cocktails are designed for use with our sample oxidizers to ensure superior performance and consistently reliable results.

Product	Product no.	Size
Carbo-Sorb E - High capacity radioactive carbo dioxide absorber	6013721	1 x 1 L
<b>MonoPhase S</b> - Highest H <sub>3</sub> counting efficiency in pure water samples up to 23%	6013109	2 x 5 L
PermaFluor E* - For counting 14CO <sub>2</sub>	6013181	1 x 1 L
samples trapped in Carbo-Sorb E	6013187	4 x 2.5 L

# Tissue solubilizers

Our tissue solubilizers are ideal for dissolving wet tissue, aqueous tissue homogenates, proteins, nucleotides, and other substances into a solution.

Product	Product no.	Size
<b>SOLVABLE™</b> - Aqueous based for wet tissue, homogenates, proteins, nucleotides	6NE9100	500 mL
<b>Soluene™ 350</b> - Strong organic-based for wet tissue, homogenates, proteins, nucleotides	6003038	500 mL
<b>Hyamine Hydroxide 10X</b> - Solubilize biological tissues and $^{14}\mathrm{CO}_2$ trapping agent	6003005	500 mL

# Autoradiography enhancers

Product	Product no.	Size
En3hance - Polyacrylamide and agarose gels, no DMSO, colorless	6NE9701	1 L
<b>Enlightning</b> - Rapid one-step enhancer for polyacrylamide gels and agarose, no DMSO	6NE9741	1 L

# Decontaminants and cleaners

Product	Product no.	Size
COUNT OFF - Liquid concentrate removing radioactive residues		1 x 2.5 L
		4 x 2.5 L
COUNT OFF Surface Cleaner - Decontaminates radioactive spills	6NE942T	6 x 22 oz pump bottles (650 mL)
<b>Pico-Kleen N</b> - Concentrated liquid cleaning and radioactive decontaminant agent	6013819	-

# Vials and caps

Our glass vials are chemically inert, making them suitable for use with aggressive reagents and solubilizers, offering good visibility and no solvent permeation. Our plastic vials deliver low background and high counting efficiency and are combustible for easier disposal.



# Miniature vials

Product	Product no.	Description
Diag Dua Wal A mal	6000252	2,000/case Economically packed. Caps packed separately
Pico Pro Vial - 4 mL	6000253	2,000/case with exclusive Anti-Static treatment
Pico Prias Vial - 6 mL	6000192	2,000/case Economically packed. Caps packed separately
Pico Prias Viai - 0 mL	6000193	2,000/case with exclusive anti-static treatment
Dies (Hang In/ Viel 4 m)	6000186	2,000/case Economically packed. Caps packed separately
Pico 'Hang-In' Vial - 6 mL	6000187	2,000/case with exclusive Anti-Static treatment
Pico Glass Vial - 7 mL	6000167	1,000/case Shrink-wrapped in 5 trays of 200 foil-lined white urea screw caps packed separately
Damie Vial 4 mal	6000592	1,000/case Shrink-wrapped in 5 trays of 200 Caps packed separately
Pony Vial - 6 mL	6000292	2,000/case Economically packed. Caps packed separately
	6000293	2,000/case with exclusive Anti-Static treatment
Polypropylene Vial for MicroBeta - 4 mL	1200-421	3,000/case Economically packed. Caps packed separately
Pico Glass Vial Caps	6000179	1,000/case White urea screw caps with foil liner

### Midsize vials

Product	Product no.	Description
11" · · · · · · · · · · · · · · · · · ·	6000480	2,000/case. Economically packed. Caps packed separately
Hinge-Cap Via - 8 mL	6000488	500/case. Shrink-wrapped in 5 trays of 100 vials each
Mid: Vial O ml	6000288	2,000/case. Economically packed. Caps packed separately
Midi-Vial - 8 mL	6000289	2,000/case. Same as above, with exclusive Anti-Static treatment
Maxi-Vial - 18 mL	6000201	1,000/case. Economically packed. Caps packed separately

# Standard vials

Product	Product no.	Description
Super-Polyethylene Vial with Glass vial thread - 20 mL	6001087	500/case Packed in 5 trays of 100 vials each with foil-lined urea screw caps on
	6008117	1,000/case Economically packed. Caps packed separately
Come Delivership and Mark	6008118	1,000/case with exclusive Anti-Static treatment
Super-Polyethylene Vial with Quick Closure - 20 mL	6000375	500/case Shrink-wrapped in 5 trays of 100 vials each. Caps packed separately
	6001075	500/case Packed in 5 trays of 100 vials each with caps on
Low-Diffusion Polyethylene Vial - 20 mL	6000477	100/case Shrink-wrapped in partitioned tray with caps on, with exclusive antistatic treatment
	6001009	500/case Packed in 5 partitioned trays of 100 vials each with caps on
	6000349	500/case Packed in 5 trays of 100 vials each with foil-lined urea screw caps on
High-Performance Glass Vial - 20 mL 6000128		500/case Shrink-wrapped in 10 trays of 50 vials each. Foil-lined urea screw caps packed separately
	6000129	500/case Shrink-wrapped in 5 trays of 100 vials. Foil-lined urea screw caps packed separately
	6000134	500/case Shrink-wrapped in 10 trays of 50 vials each. Poly-Cone lined urea crew caps packed separately

Product	Product no.	Description
Econo-Glass Vial - 20 mL	6000096	500/case Shrink-wrapped in 4 trays of 125 vials each. Foil-lined urea screw caps packed separately
Oximate Vial - 20 mL	6001095	500/case Polyethylene vials with glass vial thread. Packed in 5 trays of 100 vials each with foil-lined urea screw caps on. Special design screw for automated handling for automated handling
	6001009	Packed in 5 trays of 100 vials each with foil-lined urea screw caps on. Special design screw cap for automated handling
Glass Vial Caps	6001025	1,000/case Foil-lined urea screw caps
	6001027	1,000/case Poly-Cone lined urea screw caps
	6000249	1,000/case Poly screw caps for glass vials
	6000250	1,000/case Foil-lined urea screw caps Design for automated handling



