

# MultiTox-Fluor Multiplex Cytotoxicity Assay

The MultiTox-Fluor Multiplex Cytotoxicity Assay is a single-reagent-addition, homogeneous, fluorescent assay that measures the number of live and dead cells simultaneously in culture wells. The assay simultaneously measures cell viability and cytotoxicity by detecting two distinct protease activities. The live-cell protease activity is restricted to intact viable cells and is measured using a fluorogenic, cell-permeant peptide substrate (GF-AFC Substrate). The substrate enters intact cells where it is cleaved to generate a fluorescent signal proportional to the number of living cells. This live-cell protease activ...

[Expand to Read More »](#)



[Learn more about the Helix program](#)



Prices valid for US customers only

PRODUCT	SIZE	CONC.	CATALOG #	*LIST PRICE	QTY
MultiTox-Fluor Multiplex Cytotoxicity Assay <a href="#">Protocol</a> <a href="#">MSDS</a> <a href="#">Components</a> <a href="#">Add to Helix</a>	10ml	-	G9200	\$ 234.00	<input type="text"/> <a href="#">Add to cart</a>
MultiTox-Fluor Multiplex Cytotoxicity Assay <a href="#">Protocol</a> <a href="#">MSDS</a> <a href="#">Components</a> <a href="#">Add to Helix</a>	5 × 10ml	-	G9201	\$ 638.00	<input type="text"/> <a href="#">Add to cart</a>
MultiTox-Fluor Multiplex Cytotoxicity Assay <a href="#">Protocol</a> <a href="#">MSDS</a> <a href="#">Components</a> <a href="#">Add to Helix</a>	2 × 50ml	-	G9202	\$ 1,159.00	<input type="text"/> <a href="#">Add to cart</a>

## Storage Conditions

Store at -20°C.

For product intended use please see Patents & Disclaimers tab.

## You Also May Be Interested In



[ApoTox-Glo™ Triplex Assay](#)

[ApoLive-Glo™ Multiplex Assay](#)

[GloMax®-Multi+ Detection System with Instinct® Software](#)

[CytoTox-Fluor™ Cytotoxicity Assay](#)