



Luminescent cell viability and reporter gene assays.

Discover our Lites family of products which use luciferase for monitoring cell health or the study of transcriptional gene expression and virus life cycles. Take advantage of luminescence technology:

- Highly sensitive, wide dynamic range
- Simple and reproducible assays
- Stable at 2 - 8°C
- No DTT, eliminating toxic hazards
- Cost-effective
- Automation compatible

Illuminate your lab with a free sample

Try before you buy - request a sample to see the exceptional performance firsthand.

Qualifying products	
Part number	Product name
6016943	ATPlite 300 assay
6016736	ATPlite 1step, 10mL
6016716	neolite 10mL
6066706	twinlite 100 assay
6016796	twinlite glow 100 assay
6066726	sensilite 100 assay
6066766	britelite plus 10mL
6066756	steadylite plus 10mL
6096736	ATPlite 1glow F 10ml (Frozen) NEW
6096746	ATPlite 1glow 10ml (Lyophilized) NEW

*Terms and conditions apply



Order your free sample!

Scan or click for more information