
CERTIFICATE OF ANALYSIS

Catalog Number C014-1
Product Name NucView Red Live nucleic acid stain, 1 mM in DMSO
Appearance blue solution
Lot Number

	LOT DATA	SPECIFICATION
Absorption¹ Optical Density	0.73 at 635 nm	0.73 ± 0.05 at 635 ± 4 nm

1. Solvent: Methanol. Method: Optical Density. 200x dilution.