

CERTIFICATE OF ANALYSIS

Catalog Number	C040	
Product Name	CellView Orange CMTMR, Tetramethylrhodamine]	[5-(6)-(((4-Chloromethyl)Benzoyl)Amino)
CAS Number	323192-14-9	
Molecular Formula	C ₃₂ H ₂₈ N ₃ O ₄ Cl	
Molecular Weight	554.04	
Appearance	purple solid	
Lot Number		

	LOT DATA	SPECIFICATION
Absorption ¹ Maximum Extinction	540 nm 99,000 cm ⁻¹ M ⁻¹	540 ± 3 nm 102,000 ± 12,000 cm ⁻¹ M ⁻¹
Fluorescence ¹ Emission Maximum	565 nm	566 ± 4 nm
HPLC Purity	86 % at 254 nm	≥ 85 % at 254 nm
NMR Result	meets specification	consistent with structure

1. Solvent: Methanol.