

CERTIFICATE OF ANALYSIS

Catalog Number C022
Product Name calcein blue, AM
CAS Number 168482-84-6
Molecular Formula C₂₁H₂₃NO₁₁
Molecular Weight 465.41
Appearance off-white solid
Lot Number

	LOT DATA	SPECIFICATION
Absorption ¹		
Maximum	323 nm	322 ± 3 nm
Extinction	14,000 cm ⁻¹ M ⁻¹	13,000 ± 1,000 cm ⁻¹ M ⁻¹
Fluorescence ¹		
Emission Maximum	439 nm	437 ± 4 nm
HPLC		
Purity	95 % at 254 nm	≥ 80 % at 254 nm
NMR		
Result	meets specification	consistent with structure

1. Solvent: Methanol.