

CERTIFICATE OF ANALYSIS

Catalog Number C025
Product Name calcein, AM
CAS Number 148504-34-1
Molecular Formula C₄₆H₄₆N₂O₂₃
Molecular Weight 994.87
Appearance yellow solid
Lot Number

	LOT DATA	SPECIFICATION
Absorption ¹ Maximum Extinction	284 nm 5,600 cm ⁻¹ M ⁻¹	283 ± 3 nm 5,500 ± 500 cm ⁻¹ M ⁻¹
HPLC Purity	90 % at 254 nm	≥ 85 % at 254 nm
NMR Result	meets specification	consistent with structure

1. Solvent: Methanol.