

CERTIFICATE OF ANALYSIS

Catalog Number C026
Product Name calcein, AM 1 mg/mL solution in DMSO
CAS Number 148504-34-1
Molecular Formula $C_{46}H_{46}N_2O_{23}$
Molecular Weight 994.87
Appearance colorless solution
Lot Number

	LOT DATA	SPECIFICATION
Absorption¹ Optical Density	0.50 at 284 nm	0.50 ± 0.05 at 283 ± 3 nm

1. Solvent: Methanol. Method: Optical Density. 11× dilution.