

CERTIFICATE OF ANALYSIS

Catalog Number C052
Product Name Fluorescein phalloidin
CAS Number 84313-38-2
Molecular Formula $C_{56}H_{60}N_{10}O_{15}S_2$
Molecular Weight 1177.27
Appearance orange solid
Lot Number

	LOT DATA	SPECIFICATION
Absorption ¹ Maximum	496 nm	496 ± 3 nm
Fluorescence ¹ Emission Maximum	518 nm	518 ± 4 nm
HPLC Purity	95 % at 254 nm	≥ 90 % at 254 nm

1. Solvent: Buffer, 50 mM potassium phosphate (pH 8).