

CERTIFICATE OF ANALYSIS

Catalog Number C053
Product Name Rhodamine phalloidin
CAS Number 219920-04-4
Molecular Weight ~1250
Appearance red solid
Lot Number

	LOT DATA	SPECIFICATION
Absorption¹ Maximum	541 nm	540 ± 3 nm
Fluorescence¹ Emission Maximum	565nm	565 ± 4 nm
HPLC Purity	95 % at 254 nm	≥ 90 % at 254 nm

1. Solvent: Methanol.