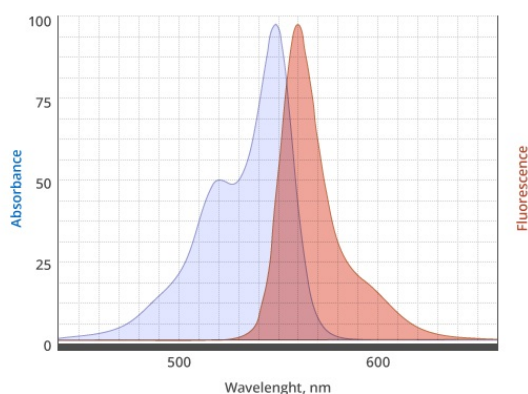


**AF555 NHS ester**

structural analog to Alexa Fluor® 555

Cat. No.	Amount
APC-003-1	1 mg
APC-003-5	5 mg



excitation and emission spectrum of AF555

For research use only!**Shipping:** shipped at ambient temperature**Storage Conditions:** store at -20 °C**Additional Storage Conditions:** store dark**Shelf Life:** 12 months after date of delivery**Molecular Weight:** 944.07 g/mol (protonated)**Purity:** > 90 % (HPLC)**Form:** solid**Solubility:** DMSO, DMF**Spectroscopic Properties:** λ_{exc} 555 nm, λ_{em} 572 nm,
 ϵ 155.0 L mmol⁻¹ cm⁻¹