

DEAC,SE [7-Diethylaminocoumarin-3-carboxylic acid, succinimidyl ester] *CAS 139346-57-9*

Catalog Number: 506

Unit Size: 25 mg

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Solid off white
Molecular Weight	358.35
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	411 nm
Emission Wavelength	472 nm

Applications

The amine-reactive 7-diethylaminocoumarin-3-carboxylic acid, succinimidyl ester can be used to create blue-fluorescent bioconjugates. It has quite strong blue fluorescence. When compared with AMCA conjugates, conjugates of the UV-light-excitable 7-dialkylaminocoumarin fluorophore have much longer-wavelength excitation and slightly longer-wavelength emission spectra (~470 nm). This fluorescent dye is quite hydrophobic, and might be used for labeling live cells.