

DEAC [7-Diethylaminocoumarin-3-carboxylic acid] *CAS 50995-74-9*

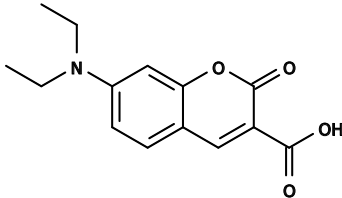
Catalog Number: 505

Unit Size: 100 mg

Product Details

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

Chemical Properties

Appearance	Solid yellow
Molecular Weight	261.27
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	411 nm
Emission Wavelength	472 nm

Applications

7-Diethylaminocoumarin-3-carboxylic acid 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).