

DDAO Galactoside

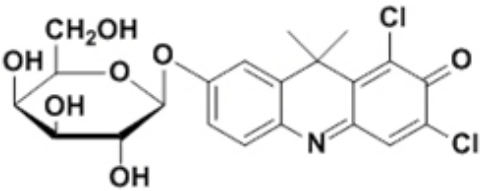
Catalog Number	Packaging Size
C281	2 mg

Storage upon receipt: -20°C, protected from light

Introduction

DDAO galactoside is a galactosidase substrate that yields a hydrolysis product that can be excited with the 633 nm laser (excitation/emission maxima ~645/660). Although the substrate itself is fluorescent (excitation/emission maxima ~460/610 nm), the difference between the substrate's and the hydrolysis product is over 200 nm, along the two species to be easily distinguished.

Specifications

Label:	DDAO	
Ex/Em:	646/659 nm	
Detection Method:	Fluorescent	
Molecular Formula:	C ₂₁ H ₂₁ Cl ₂ NO ₇	
Molecular Weight:	470.30	
CAS Number:	503178-95-8	
Storage Conditions:	-20°C, protected from light	
Shipping Condition:	Room Temperature	

Applications

Galactosidase Substrate

References:

1. [In vivo imaging of beta-galactosidase activity using far red fluorescent switch.](#)
Tung CH, Zeng Q, Shah K, Kim DE, Schellingerhout D, Weissleder R
Cancer Res (2004) 64:1579-1583
2. [Enzymatic detection of protein translocation.](#)
Wehrman TS, Casipit CL, Gewertz NM, Blau HM
Nat Methods (2005) 2:521-527
3. [Oxygen sensitivity of reporter genes: implications for preclinical imaging of tumor hypoxia.](#)
Cecic I, Chan DA, Sutphin PD, Ray P, Gambhir SS, Giaccia AJ, Graves EE
Mol Imaging (2007) 6:219-228