

DDAO phosphate

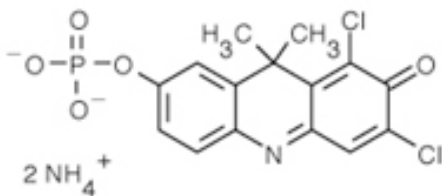
Catalog Number	Packaging Size
C276	2 mg

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

Introduction

The phosphatase substrate **DDAO phosphate** [9H-(1,3-dichloro-9,9-dimethylacridin-2-one-7-yl) phosphate, diammonium salt] yields a hydrolysis product that can be excited with the 633 nm laser (excitation/emission maxima ~646/659). Although the substrate itself is fluorescent (excitation/emission maxima ~460/610 nm), the difference between the substrate's excitation maximum and that of its hydrolysis product is over 200 nm, allowing the two species to be easily distinguished.

Specifications

Label:	DDAO	
Ex/Em:	646/659 nm	
Detection Method:	Fluorescent	
Molecular Formula:	C ₁₅ H ₁₈ Cl ₂ N ₃ O ₅ P	
Molecular Weight:	422.20	
CAS Number:	500883-59-0	
Storage Conditions:	-20°C, protect from light	
Shipping Condition:	Room Temperature	

Applications

Phosphatase Substrate

References:

1. Simultaneous trichromatic fluorescence detection of proteins on Western blots using an amine-reactive dye in combination with alkaline phosphatase- and horseradish peroxidase-antibody conjugates. Simultaneous trichromatic fluorescence detection of proteins on Western blots using an amine-reactive dye in combination with alkaline phosphatase- and horseradish peroxidase-antibody conjugates. Martin K, Hart C, Liu J, Leung WY, Patton WF. Proteomics (2003) 3:1215-1227
2. Single-molecule detection technologies in miniaturized high-throughput screening: fluorescence intensity distribution analysis. Single-molecule detection technologies in

miniaturized high-throughput screening: fluorescence intensity distribution analysis.
Haupts U, Rudiger M, Ashman S, Turconi S, Bingham R, Wharton C, Hutchinson J,
Carey C, Moore KJ, Pope AJ
J Biomol Screen (2003) 0: 19-19

3. Global quantitative phosphoprotein analysis using multiplexed proteomics
technology. Global quantitative phosphoprotein analysis using multiplexed proteomics
technology.
Steinberg TH, Agnew BJ, Gee KR, Leung W-Y, Goodman T, Schulenberg B, Hendrickson
J, Beechem JM, Haugland RP, Patton WF
Proteomics (2003) 3: 1128-1128