

# DDAO Phosphate (1 mg)

DDAO-phosphate is a long-wavelength, phosphatase substrate for both solution-based assays and the amplified detection of specific targets on blots. Hydrolysis of the phosphate bond yields a brightly red-fluorescent product with absorption/emission maxima of ~646/659 nm that is efficiently excited by the 633 nm spectral line of the He–Ne laser. The shift of emission maxima between DDAO-phosphate and the dephosphorylated product is larger than 200 nm and permits both species to be easily distinguished. DDAO phosphate has good water solubility, a low  $K_M$  and a high turnover rate.

## Ordering Information

[Order Online »](#)

JBS-APC-001-1	1mg
---------------	-----

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 9H-(1,3-Dichloro-9,9-dimethylacridin-2-one-7-yl) phosphate, Bis(triethylammonium) salt

**Appearance** solid

**CAS** 500883-59-0

**Emission Maximum** 628 nm

**Excitation Maximum** 477 nm

**Formula**  $C_{15}H_{12}Cl_2NO_5P$  (free acid)

**MW** 388.14 (free acid)

**Purity** > 95 % (HPLC)

**Technical Info / Product Notes** For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)