

DMEQ-TAD

DMEQ-TAD is a reagent for fluorescence-labeling of Vitamin D and its use in assaying Vitamin D metabolites.

Citations: 1

[View Online »](#)

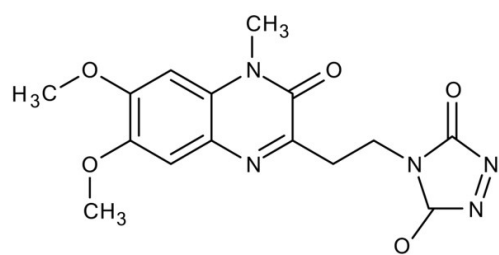
Ordering Information

[Order Online »](#)

ENZ-CHM249-0100	100mg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Store under nitrogen in the dark.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	4-[2-(6,7-Dimethoxy-4-methyl-3-oxo-3,4-dihydroquinoxaliny)ethyl]-1,2,4-triazoline-3,5-dione
Appearance	Bright pink solid.
CAS	132788-52-4
Formula	$C_{15}H_{15}N_5O_5$
Identity	Determined by NMR.
MW	345.3
Purity	≥98% (TLC)
Solubility	Soluble in dichloromethane, acetone or chloroform.

Last modified: May 29, 2024



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-
fr@enzolifesciences.com](mailto:info-fr@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)