

Digoxigenin NHS ester

Activated ester that reacts with amino groups

Digoxigenin is a non-radioactive immune-tag, widely used for *in situ* hybridization and protein labeling. Pairing with anti-digoxigenin antibodies, digoxigenin tag allows the detection of interested biomolecules with high affinity and specificity. Digoxigenin NHS ester reacts with amino groups within biomolecules (nucleic acid, proteins, etc) and enables the attachment of digoxigenin into these molecules.

Citations: 4

[View Online »](#)

Ordering Information

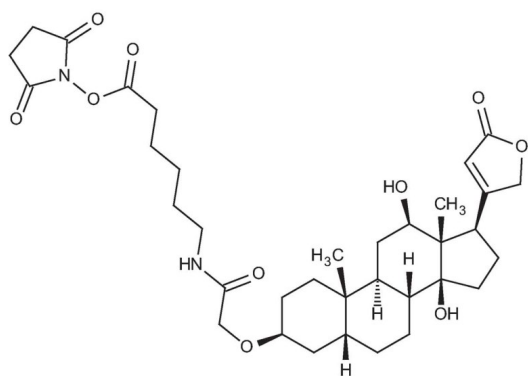
[Order Online »](#)

ENZ-45022-0001	1mg
ENZ-45022-0005	5mg

Manuals, SDS & CofA

[View Online »](#)

- Used for labeling of proteins and oligonucleotides
- Allows to immuno-tag molecules for detection using anti-digoxigenin antibodies with high affinities and specificity



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Handling	Store in a dry place.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Off-white solid.
CAS	129273-26-3
Formula	$C_{35}H_{50}N_2O_{10}$
MW	658.78
Purity	≥80% (HPLC)
Solubility	Soluble in DMSO or DMF.



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)