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

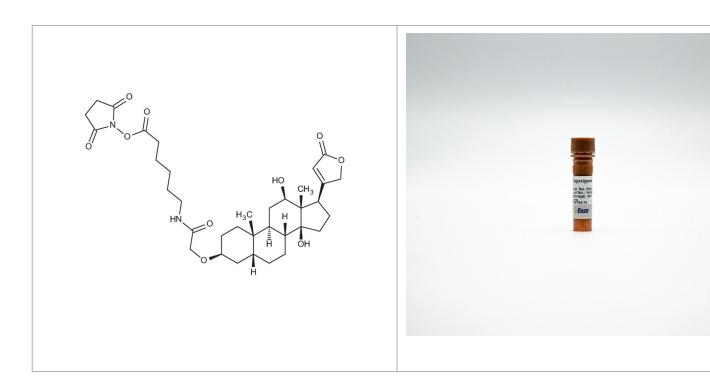
Order Online »

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 antidigoxigenin 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.

Last modified: May 29, 2024

