

Digoxigenin C2 acid

Catalog Number: 3509

Unit Size: 5 mg

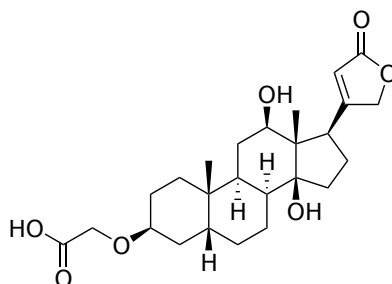
Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	24 months upon receiving

Chemical Properties

Appearance	Solid
Molecular Weight	448.56
Soluble In	DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

Digoxigenin (DIG) is a steroid found exclusively in the flowers and leaves of the plants *Digitalis purpurea*, *Digitalis orientalis*, and *Digitalis lanata* (foxgloves). It is attached to sugars, forming glycosides such as digoxin and lanatoside C. In biological detections, DIG serves as a hapten, a small molecule with high antigenicity. Similar to other popular haptens like 2,4-dinitrophenol (DNP), biotin, and fluorescein, DIG can be readily conjugated into biomolecules (such as proteins and nucleic acids) for a variety of biological detections and tests. It is useful for immunohistochemistry and in situ hybridization. For example, DIG can be conjugated to a single species of RNA nucleoside triphosphate (typically uridine) during RNA synthesis, creating sensitive non-radioactive riboprobes to detect nucleic acids in plants and animals. These riboprobes can detect as little as 1 µg of plasmid DNA. Anti-digoxigenin antibodies have high affinities and are specifically used in various biological immuno-assays, including ELISA.