

# Genistin

## TdT inhibitor

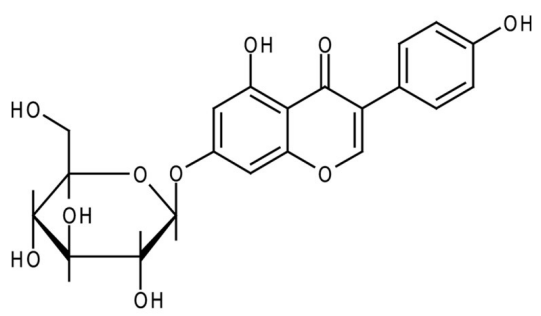
Glucoside of genistein (Prod. No. ALX-350-006) found in soy beans. Useful as a negative control for genistein and other tyrosine kinase inhibitors. Selective inhibitor of terminal deoxyribonucleotidyltransferase (TdT). Displays antioxidant and anticarcinogenic properties.

Citations: 7

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** -20°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Genistein-7-O-glucoside

**Appearance** Off-white to light yellow powder.

**CAS** 529-59-9

**Couple Target** TdT

**Couple Type** Inhibitor

**Formula**  $C_{21}H_{20}O_{10}$

**MI** **14:** 4391

**MW** 432.4

**Purity** ≥98% (HPLC, TLC)

**RTECS** DJ3093000

**Solubility** Soluble in DMSO or dilute aqueous base; insoluble in water.

**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-  
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)