

IPTG (animal free) (free of dioxane)

β -galactosidase activator

IPTG mimicks allolactose, a metabolite of lactose. IPTG is used to trigger protein expression where the gene is under the control of the *lac* operon. IPTG can be used along X-Gal (Prod. No. ALX-582-002) to detect the activity of the *lac* gene in cloning experiments.

Citations: 1

[View Online »](#)

Ordering Information

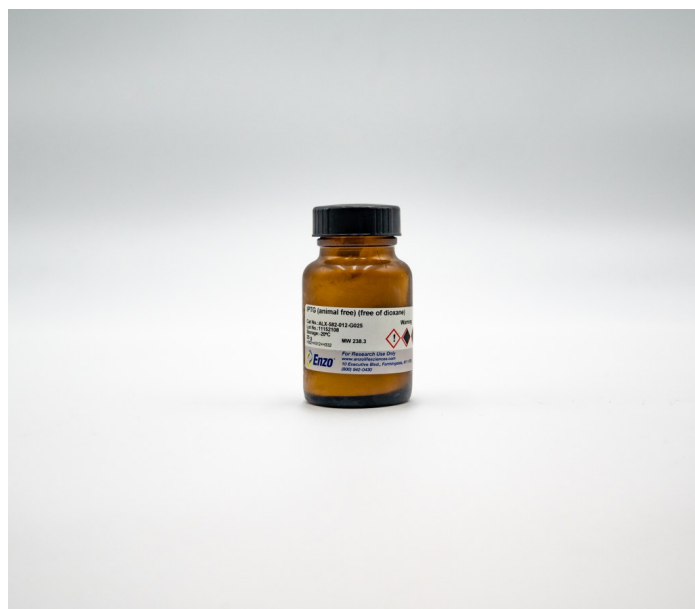
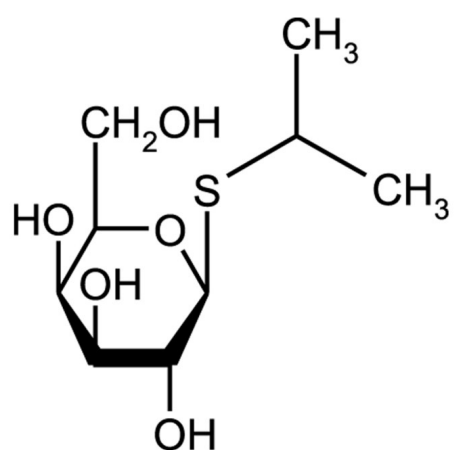
[Order Online »](#)

ALX-582-012-G005	5g
ALX-582-012-G025	25g

Manuals, SDS & CofA

[View Online »](#)

- Inducer of β -Galactosidase
- Used in the stimulation of β -Galactosidase in cellular systems in which dioxane would disrupt normal cell function
- Contains no animal-derived products



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	1-methylethyl 1-thio-β-D-galactopyranoside, Isopropyl β-D-thiogalactoside
Appearance	White to off-white powder.
CAS	367-93-1
Couple Target	Galactosidase
Couple Type	Activator
Formula	$C_9H_{18}O_5S$
Identity	Identity determined by IR.
MW	238.3
Purity	≥98% (HPLC)
Solubility	Soluble in 100% methanol or water (5%).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com