

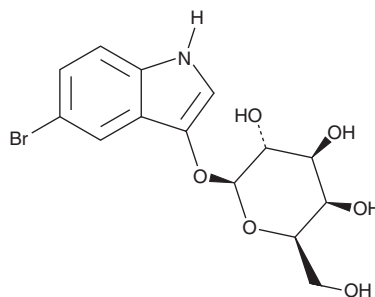
PRODUCT INFORMATION



5-bromo-3-indolyl-β-D-Galactopyranoside

Item No. 35299

CAS Registry No.: 97753-82-7
Formal Name: 5-bromo-1H-indol-3-yl β-D-galactopyranoside
Synonym: Bluo-Gal
MF: C₁₄H₁₆BrNO₆
FW: 374.2
Purity: ≥98%
UV/Vis.: λ_{max}: 228 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

5-bromo-3-indolyl-β-D-Galactopyranoside (Bluo-Gal) is supplied as a solid. A stock solution may be made by dissolving the Bluo-Gal in the solvent of choice, which should be purged with an inert gas. Bluo-Gal is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of Bluo-Gal in these solvents is approximately 2 and 1 mg/ml, respectively.

Description

Bluo-Gal is a cell-permeable chromogenic substrate for β-galactosidase.¹ Upon enzymatic hydrolysis, it generates an indigo-colored precipitate that can be used to measure β-galactosidase activity. Bluo-Gal has been used to detect the expression of β-galactosidase in muscle from transgenic mice expressing *E. coli lacZ*, the gene that encodes β-galactosidase.

Reference

1. Gioglio, L., de Angelis, M.G.C., Boratto, R., *et al.* An improved method for β-galactosidase activity detection on muscle tissue. A light and electron microscopic study. *Ann. Anat.* **184**(2), 153-157 (2002).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/18/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM