

Glowing products for science

IPTG

IPTG (Isopropylthio- β -galactoside) is an inducer of β -galactosidase activity in bacteria and is suitable for use with X-gal or bluo-gal to detect lac gene activity in cloning procedures.



Product attributes

Call us: 800-304-5357

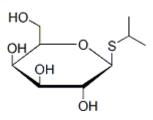
CAS number

367-93-1

Product Description

IPTG (Isopropylthio- β -galactoside) is an inducer of β -galactosidase activity in bacteria and is suitable for use with X-gal or bluo-gal to detect lac gene activity in cloning procedures.

- White solid soluble in DMSO
- Store desiccated at 4°C and protect from light
- C₉H₁₈O₅S
- MW: 238.3



This datasheet was generated on December 2, 2025 at 10:31:06 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/iptg/