

IPTG

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



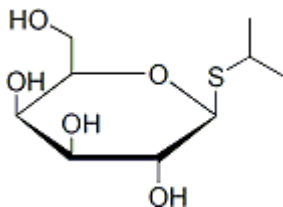
Product attributes

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 blue-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 June 20, 2026 at 07:25:38 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/iptg/>