



Isopropyl β -D-thiogalactopyranoside (IPTG)
 β -galactosidase Inducer
Molecular Weight 238.31

Catalog Number 170-3201

Contents	1 gram
Storage	-20 °C
Use	IPTG is a gratuitous inducer of bacterial β -galactosidase. This product has been specially qualified for the detection of phage plaques or plasmid/phagemid containing bacterial colonies expressing protein fused to β -galactosidase.
Quality assurance	IPTG has been tested functionally with <i>E. coli</i> MV1190 infected with M13 phage. Blue plaques form only in the presence of IPTG and X-gal (5-bromo-4-chloro-3-indolyl- β -D-galactopyranoside), a substrate of β -galactosidase.

This product is intended for research use only.