

## PRODUCT INFORMATION

# X-Gal Solution, ready-to-use

5-bromo-4-chloro-3-indolyl- $\beta$ -D-galactopyranoside

#\_ \_ \_

Lot: \_ \_ \_ Expiry Date: \_ \_

Formula:  $C_{14}H_{15}BrClNO_6$   
Molecular Weight: 408.6  
Concentration: 20 mg/ml

**Store at 4°C (or at room temperature)  
in the original vial.**

## Description

X-Gal is a noninducing chromogenic substrate of  $\beta$ -galactosidase, which hydrolyzes X-Gal forming an intense blue precipitate. X-Gal is most frequently used in conjunction with IPTG in blue/white colony screening to detect recombinants (white) from non-recombinants (blue).

X-Gal Solution, ready-to-use, is a stabilized non-toxic 20 mg/ml solution of X-Gal.

## Application

Direct use in blue/white colony screening experiments.

## CERTIFICATE OF ANALYSIS

Tested in the blue/white cloning assay.

Quality authorized by:  Jurgita Zilinskiene

## RECOMMENDATIONS FOR USE

### For individual LB plates

- Let LB plates warm at room temperature under UV.
- Spread out evenly 40 µl of the X-Gal Solution, ready-to-use on the Petri dish surface.

### For batch use

- Add directly 2 µl of X-Gal Solution, ready-to-use, to every milliliter of molten LB agar (temperature 50°C or less).
- Mix well and pour.
- Let cool.

### Note

The appearance of pale blue or light yellow color of X-Gal Solution, ready-to-use, does not have influence on its quality.

## SAFETY INFORMATION



### X-Gal Solution, ready-to-use

**Xi** Irritant

### Risk phrases

R10 Flammable.

R36/38 Irritating to eyes and skin.

### Safety phrases

S7 Keep container tightly closed.

S16 Keep away from sources of ignition - No smoking.

S24/25 Avoid contact with skin and eyes.

S26 In case of contact with eyes, rinse immediately with plenty of water and seek medical advice.

S37 Wear suitable gloves.

S60 This material and its container must be disposed of as hazardous waste.

### PRODUCT USE LIMITATION

This product is developed, designed and sold exclusively *for research purposes and in vitro use only*. The product was not tested for use in diagnostics or for drug development, nor is it suitable for administration to humans or animals. Please refer to [www.thermoscientific.com/onebio](http://www.thermoscientific.com/onebio) for Material Safety Data Sheet of the product.

© 2012 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.