

46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352

www.biotium.com

PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Rose-β-D-Gal (6-chloro-3-indoxyl-β-D-galactopyranoside)

10013 **CATALOG NUMBER:**

MOLECULAR C₁₄H₁₆ClNO₆ **INFORMATION:** Mwt: 329.74.

[138182-21-5]

PROPERTIES:

Color & Form White solid. **Purity** \geq 98% by TLC **Solubility** Soluble in MeOH

STORAGE AND HANDLING:

Store desiccated at ≤ 4 °C upon receipt. Protect from light, especially in solution.

APPLICATION: Chromogenic substrate for β -galactosidase. The product is similar to X-gal, but

generates an intense pink colored ($\lambda_{max} = 542 \text{ nm}$) precipitate on enzymatic hydrolysis. Protocol for using the material is similar to that for X-gal (See references listed below).

Ref: 1) Molecular Cloning: A laboratory Manual, Cold Spring Harbor Laboratory, B. 14, p186(1989);

2) Biotechniques, 27, 438(1999);

3) Oncogen, 10, 2323(1995);

4) Devel. Biol., 161, 77(1994);

5) J. Histochem. Cytochem., 42, 1299(1994);

6) BioTechniques, 15, 974(1993);

7) BioTechniques, 15, 292(1993).

TOXICITY: Unknown

Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on FIRST AID:

clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing,

move individual to fresh air and seek medical advice immediately.

Disclaimer: Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.