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:** X-Gal (5-Bromo-4-chloro-3-indoxyl-beta-D-galactopyranoside)

CATALOG #: 10011/10011-1/10011-2

**MOLECULAR**  $C_{14}H_{15}BrCINO_6$  **INFORMATION:** MW: 408.61 [7240-90-6]

OH OH OH OH NH

**PROPERTIES:** 

Color & Form White, crystalline power

**Purity**  $\geq$  99% by HPLC

**Soluble** in DMSO or DMF

STORAGE AND HANDLING:

Store desiccated at -20°C. Hygroscopic! Keep tightly closed.

**APPLICATION:** X-gal is the most widely used chromogenic substrate for  $\beta$ -galactosidase. Enzyme

hydrolysis of the substrate yields a dark blue colored precipitate ( $\lambda_{abs} \sim 615 \text{ nm}$ )

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.

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

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.