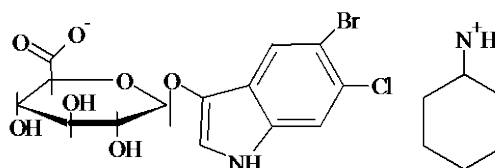


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: **Red-β-D-GlcU** (Magenta-β-D-GlcUA·CHX, 5-bromo-6-chloro-3-indolyl-β-D-glucuronide, cyclohexylammonium salt)

CATALOG NUMBER: 10018

MOLECULAR INFORMATION: C₂₀H₂₆BrClN₂O₇
Mwt: 521.8
[144110-43-0]

**PROPERTIES:**

Color & Form White solid
Purity ≥ 98% by TLC
Solubility Soluble in DMF or DMSO

STORAGE AND HANDLING:

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

APPLICATION:

This product is similar to X-GlcU, but the product of the substrate is a red precipitate, which is easier to detect against the background of plant cells.

Ref: *BioTechniques* 19(3), 352 (1995)

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.