

PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Chromogenic β -Glucuronidase Substrate Sampler Kit

CATALOG #: 10025

MOLECULAR INFORMATION: See Individual Components

PROPERTIES:

Color & Form See Individual Components

Purity See Individual Components

Solubility See Individual Components

STORAGE AND HANDLING: See Individual Components

APPLICATION: This kit contains 5 mg each of the following three chromogenic β -glucuronidase substrates: X-GlcU, CHA (#10016), Red- β -D-GlcU, CHA (#10018), and Rose- β -D-GlcU, CHA (#10019)

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: <i>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.</i>
--