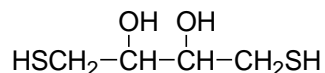


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: DTT (dithiothreitol)**CATALOG #** 91050**MOLECULAR INFORMATION:** C₄H₁₀O₂S₂
MWt: 154.2
[3483-12-3]**PROPERTIES:****Color & Form** White solid
Solubility Soluble in water (5% solution in water)
Absorption/Emission Extinction Coefficient**STORAGE AND HANDLING:** Store at ≤ -20°C**APPLICATION:** DTT is a reducing agent commonly used to reduce cystine disulfides to free thiols.Ref.: 1) *Anal Biochem* **273**, 73(1999); 2) *Methods Enzymol* **143**, 246(1987)**TOXICITY:** Not established. Not listed by NTP, IARC or OSHA.

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>
--