CTC (5-Cyano-2,3-ditolyl tetrazolium chloride)

5-Cyano-2,3-ditolyl tetrazolium chloride (CTC) has been used to evaluate the respiratory activity of many bacterial populations as well as populations undergoing drug efficacy evaluations. Healthy cells respiring via the electron transport chain will absorb and reduce CTC into an insoluble, red fluorescent formazan product.

I HIND I HIND I COLD I

Product attributes

CAS number 102568-47-8

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

5-Cyano-2,3-ditolyl tetrazolium chloride (CTC) has been used to evaluate the respiratory activity of many bacterial populations as well as populations undergoing drug efficacy evaluations. Healthy cells respiring via the electron transport chain will absorb and reduce CTC into an insoluble, red fluorescent formazan product. Cells not respiring or respiring at slower rates will reduce less CTC, and consequently produce less fluorescent product, giving a semi-quantitative estimate of healthy vs. unhealthy bacteria.

- Pale yellow solid soluble in water
- C₁₆H₁₄CIN₅
- MW: 311.8
- [102568-47-8]

This datasheet was generated on November 12, 2025 at 09:19:55 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/ctc-5-cyano-23-ditolyl-tetrazolium-chloride/