

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



Product attributes

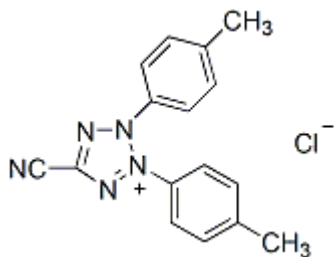
CAS number	102568-47-8
------------	-------------

5-Cyano-2,3-ditoly 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.

Product Description

5-Cyano-2,3-ditoly 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
- $C_{16}H_{14}ClN_5$
- MW: 311.8
- [102568-47-8]



This datasheet was generated on June 10, 2026 at 03:24:19 AM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/ctc-5-cyano-23-ditoly-tetrazolium-chloride/>