

BisEA

BisEA (Bis-(2-(Methanethiosulfonato)ethyl)amine, hydrochloride) is a positively charged MTS cross-linker.

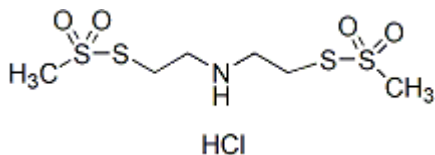


Product attributes

Product Description

BisEA (Bis-(2-(Methanethiosulfonato)ethyl)amine, hydrochloride) is a positively charged MTS cross-linker.

- Off-white solid soluble in DMF or DMSO
- Store desiccated at -20 °C
- $C_6H_{16}ClNO_4S_4$
- MW: 329.91



References

1. PNAS 98(18), 10055 (2001).

This datasheet was generated on June 15, 2026 at 01:05:48 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/bisea-bis-2-methanethiosulfonatoethylamine-hydrochloride/>