

AEAETS

AEAETS (2-Aminoethyl 2-aminoethanesulfonate dihydrochloride salt) is a doubly positively charged alkanethiosulfonate.

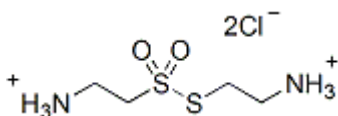


Product attributes

Product Description

AEAETS (2-Aminoethyl 2-aminoethanesulfonate dihydrochloride salt) is a doubly positively charged alkanethiosulfonate.

- White solid soluble in water, DMSO and DMF
- Store at -20 °C
- C₄H₁₄Cl₂N₂O₂S₂
- MW: 257



References

1. J Gen Physiol 115, 93 (2000).
2. Meth Enzymol 293, 123 (1998).

This datasheet was generated on June 3, 2026 at 11:32:27 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/aeaets-2-aminoethyl-2-aminoethanesulfonate-dihydrochloride-salt/>