

EDC (EDAC)

EDC (or EDAC) (1-(3-Dimethylaminopropyl-3-ethylcarbodiimide, hydrochloride)) has recently been found to be very useful for fixing *in situ* chelators, including the fluorescent ion indicators for histological studies following physiological experiments. EDC is also a widely used reagent to activate carboxy groups for amine coupling.



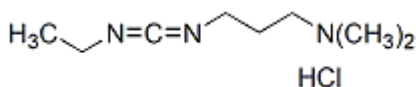
Product attributes

CAS number	25952-53-8
------------	------------

Product Description

EDC (or EDAC) (1-(3-Dimethylaminopropyl-3-ethylcarbodiimide, hydrochloride)) has recently been found to be very useful for fixing *in situ* chelators, including the fluorescent ion indicators for histological studies following physiological experiments. EDC is also a widely used reagent to activate carboxy groups for amine coupling.

- White solid soluble in water
- Store desiccated at -20 °C
- C₈H₁₈ClN₃
- MW: 191.7
- [25952-53-8]



References

1. Cell Calcium 21(3),175 (1997).
2. Tetrahedron Lett 34, 7685 (1993).

This datasheet was generated on January 22, 2026 at 11:54:45 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/edc-edac-1-3-dimethylaminopropyl-3-ethylcarbodiimide-hydrochloride/>