

Probenecid, Sodium Salt, Water Soluble

Probenecid is commonly used to inhibit organic-anion transporters located in the cell membrane. Such transporters can extrude dyes and indicators and thus contribute to low signal or high background in assays due to poor cellular retention of dyes.



Product attributes

CAS number	23795-03-1
------------	------------

Product Description

Probenecid is commonly used to inhibit organic-anion transporters located in the cell membrane. Such transporters can extrude dyes and indicators and thus contribute to low signal or high background in assays due to poor cellular retention of dyes. The use of probenecid to block the efflux of intracellular dyes was first described by Di Virgilio et al. (1990), and it has been used with a wide range of anionic dyes and conjugates. The commonly used free acid form of probenecid is difficult to dissolve, requiring 1M NaOH to prepare a solution. Our water-soluble probenecid sodium salt dissolves quickly in assay buffer and eliminates the need for handling caustic NaOH. Each vial is directly dissolved in 1 mL buffer to make a 250 mM stock solution. Typical working concentrations in cell-based assays are 1-2.5 mM.

- Colorless solid soluble in water (77 mg/mL of water)
- $C_{13}H_{18}NNaO_4S$
- MWt: 307.34
- Store desiccated at room temperature
- [23795-03-1]

References

1. PLoS ONE (2014), doi:10.1371/journal.pone.0110772

This datasheet was generated on January 19, 2026 at 03:19:23 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/probenecid-sodium-salt-water-soluble/>