

46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352

www.biotium.com Revised: April 4, 2012

PRODUCT INFORMATION

PRODUCT NAME: Probenecid, sodium salt *water soluble*

CATALOG #: 50027

SIZE: 10 x 77 mg

MOLECULARC13H18NO4SNaINFORMATION:MWt: 307

PROPERTIES:

- Color & Form White solid
- Solubility 77 mg/ml water

INTRODUCTION: Probenecid is commonly used to inhibit organic-anion transporters located in the plasma membrane. Such transporters can extrude dyes and indicators and thus contribute to poor loading or high background signal in assays that depend on the intracellular retention of dyes or indicators. The use of probenecid to block the efflux of intracellular dyes was first described by Di Virgilio et al. (1990), and subsequently it has been used with a wide range of anionic dyes and conjugates. The commonly used free acid form of probenecid has poor aqueous solubility and requires the addition of 1M NaOH to dissolve in water. Our water soluble probenecid dissolves readily in water and eliminates the need to handle caustic NaOH.

STORAGE ANDStore dry probenecid desiccated at room temperature. Product is stable for at least one
year from date of receipt when stored as recommended.

Dissolve the contents of one vial in 1 mL of water or buffer to obtain a 250 mM stock solution. Stock solutions can be stored at -20°C for up to 6 months. Typical working concentrations in cell-based assays range between 1-2.5 mM.

Ref: Di Virgilio F., et.al Cell Calcium, 1990, 11, 57.

Disclaimer: Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.