

PRODUCT DATA SHEET

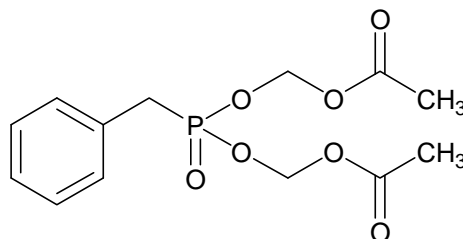


PRODUCT: Benzylphosphonic acid-(AM)₂
Benzylphosphonic acid bis acetoxymethyl ester

CATALOG NO.: EI-245

LOT NO.: temp

STRUCTURE:



PHYSICAL APPEARANCE: White solid (m.p. 38-39°C)

MOLECULAR FORMULA: C₁₃H₁₇O₇P

MOLECULAR WEIGHT: 316.3

PURITY: 98%(TLC: 5% MeOH/methylene chloride; R_f = 0.52)

SOLUBILITY: Soluble in ethanol (20 mg/mL) and DMSO (20 mg/mL)

STORAGE: Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.

APPLICATION NOTES: Avoid alkaline media (pH>8). Cell permeable analog of benzylphosphonic acid. Converted to the active phosphonic acid via intracellular esterase-catalyzed cleavage of the AM esters.

The pharmacological and toxicological properties of this product have not been fully investigated. Exercise caution in use and handling. This product must not be used in humans.