

PRODUCT DATA SHEET



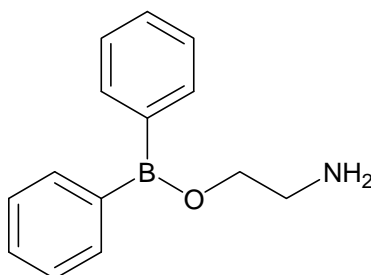
PRODUCT: 2-APB
2-Aminoethoxydiphenylborate

CAS NO.: 524-95-8

CATALOG NO.: CA-238

LOT NO.:

STRUCTURE:



PHYSICAL APPEARANCE: White solid

MOLECULAR FORMULA: $C_{14}H_{16}BNO$

MOLECULAR WEIGHT: 225.1

PURITY: 98% (TLC: 2% Methanol/methylene chloride; $R_f = 0.60$)

SOLUBILITY: Soluble in: DMSO (>25 mg/mL); ethanol (>25 mg/mL)

STORAGE: Store as supplied, at room temperature for up to 1 year. Store solutions at $-20^{\circ}C$ for up to 3 months.

APPLICATION NOTES: 2-APB activates TRPV1, TRPV2 and TRPV3 but not TRPV4, TRPV5 and TRPV6¹. It blocks TRPC5 ($IC_{50}=20 \mu M$)², TRPC6¹, TRPM3² and TRPM8¹. It potentiates the effect of capsaicin, protons and heat on TRPV1¹. It was originally introduced as an IP₃ receptor blocker and subsequently as an inhibitor of store-operated Ca^{2+} entry ($75 \mu M$)³.

REFERENCES: 1. H.-Z. Hu et al. *J. Biol. Chem.* 2004 **279** 35741
2. S.-Z. Xu et al. *Br. J. Pharmacol.* 2005 **145** 405
3. M.D. Bootman et al. *FASEB J.* 2002 **16** 1145

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.