

PRODUCT DATA SHEET



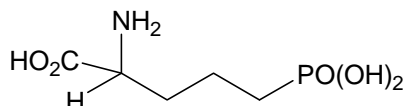
PRODUCT: DL-2-Amino-5-phosphonovaleric acid
AP5

CAS NO: 76326-31-3

CATALOG NO.: CA-400

LOT NO.: temp

STRUCTURE:



PHYSICAL APPEARANCE: White solid (m.p. 246-248°C)

MOLECULAR FORMULA: C₅H₁₂NO₅P

MOLECULAR WEIGHT: 197.1

PURITY: 99%

SOLUBILITY: Soluble in water

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

APPLICATION NOTES: Potent NMDA antagonist.^{1,2}

REFERENCES: 1. J.C.Watkins *et al. Ann.Rev.Pharmacol.Toxicol.* 1981 **21** 165
2. J.J.Hansen and P.Krogsgaard-Larsen *Med.Res.Rev.* 1990 **10** 55

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.