

PRODUCT DATA SHEET

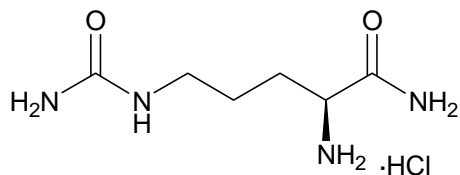


PRODUCT: FrPbAll
L-2-Amino-5-ureidopentanamide

CATALOG NO.: EA-150

LOT NO.: A11006

STRUCTURE:



PHYSICAL APPEARANCE: Off-white solid

MOLECULAR FORMULA: $C_6H_{14}N_4O_2 \cdot HCl$

MOLECULAR WEIGHT: 210.7

PURITY: 98% (NMR)

SOLUBILITY: Soluble in: Water (>25 mg/mL); DMSO (10 mg/mL); Ethanol (25 mg/mL)

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

APPLICATION NOTES: FrPbAll is a component of the *Parawixia bistriata* spider venom. It has been shown to be a selective inhibitor of GABA and glycine uptake ($IC_{50}=65 \mu M$ and $62 \mu M$ respectively) with little or no effect on monoamine or glutamate transporters¹. It produces neuroprotective effects in a rat retinal ischemia/reperfusion model displaying remarkable reduction in cell death¹. It also abolishes convulsive tonic-clonic seizures induced by bicuculline (BIOMOL Cat. # EA-149) in rats². Active *in vivo*. Crosses the blood-retina barrier.

REFERENCES: 1. R.O. Beleboni *et al. Mol. Pharmacol.* 2006 **69** 1998
2. M.A.R. Cairrao *et al. Pharm. Biol.* 2002 **40** 472

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.