

Alkaline Phosphatase

Product Number:	11900
Product Name:	Alkaline Phosphatase
E.C. Number:	3.1.3.1
CAS Number:	90031-78-9
EINECS Number:	232-631-4
Systematic Name:	Orthophosphoric-monoester phosphohydrolase (alkaline optimum)
Source:	Calf Intestine
Form:	In 50% glycerol containing PBS, 1 mM MgCl ₂ and 0.1 mM ZnCl ₂ , pH 7.6
Solubility:	Soluble in distilled water or dilute buffer
Activity:	5,000U
Unit Definition:	The amount of enzyme which liberates one micromole of p-nitrophenol from p-nitrophenyl-phosphate per minute at 37°C, pH 9.8 in diethanolamine buffer.
Storage Conditions:	2-8°C. DO NOT FREEZE