INSTRUCTIONS

1-Step PNPP



37621		0755.2
Number	Description	
37621	1-Step PNPP, 100mL, contains a non-mercury, non-azide preservative	
	Storage: Upon receipt store product at 4°C. Product is shipped at ambient temperature.	

Introduction

Thermo Scientific 1-Step PNPP (*p*-nitrophenyl phosphate disodium salt) is a widely used substrate for detecting alkaline phosphatase in ELISA-based applications. The reaction produces a water-soluble yellow product that absorbs light at 405nm. 1-Step PNPP is supplied pre-mixed with substrate buffer and is ready to use.

Procedure for using 1-Step PNPP

- 1. Equilibrate the 1-Step PNPP to room temperature.
- 2. Gently mix the 1-Step PNPP end over end.
- 3. Add 100µL of the 1-Step PNPP to each well. Mix thoroughly by gently agitating the plate.
- 4. Incubate plate at room temperature for 15-30 minutes or until sufficient color develops.
- 5. Add 50µL of 2N NaOH to stop the reaction. Mix thoroughly by gently agitating the plate.
- 6. Measure absorbance of each well at 405nm.

Related Thermo Scientific Products

31493	EZ-Link [™] Maleimide Activated Alkaline Phosphatase Kit
21323	Streptavidin, Alkaline Phosphatase Conjugated, 3mg
15075	Reagent Reservoirs, 200/pkg.
15082	Microtube Racked System, 960 tubes
15036	Sealing Tape for 96-Well Plates, 100/pkg.
37538	StartingBlock [™] (PBS) Blocking Buffer, 1L
37542	StartingBlock (TBS) Blocking Buffer, 1L

General References

- Feller, A.C., *et al.* (1983). Combined immunohistochemical staining for surface IgD and T-lymphocyte subsets with monoclonal antibodies in human tonsils. *Histochem J* 15:557-62.
- Yolken, R.H. (1982). Enzyme immunoassays for the detection of infectious agents in body fluids: Current limitations and future prospects. *Rev Infect Dis* **4:**35-68.
- Malik, N.J. and Daymon, M.E. (1982). Improved double immunoenzyme labeling using alkaline phosphatase and horseradish peroxidase. *J Clin Path* **35**:1092-4.



This product ("Product") is warranted to operate or perform substantially in conformance with published Product specifications in effect at the time of sale, as set forth in the Product documentation, specifications and/or accompanying package inserts ("Documentation") and to be free from defects in material and workmanship. Unless otherwise expressly authorized in writing, Products are supplied for research use only. No claim of suitability for use in applications regulated by FDA is made. The warranty provided herein is valid only when used by properly trained individuals. Unless otherwise stated in the Documentation, this warranty is limited to one year from date of shipment when the Product is subjected to normal, proper and intended usage. This warranty does not extend to anyone other than the original purchaser of the Product ("Buyer").

No other warranties, express or implied, are granted, including without limitation, implied warranties of merchantability, fitness for any particular purpose, or non infringement. Buyer's exclusive remedy for non-conforming Products during the warranty period is limited to replacement of or refund for the non-conforming Product(s).

There is no obligation to replace Products as the result of (i) accident, disaster or event of force majeure, (ii) misuse, fault or negligence of or by Buyer, (iii) use of the Products in a manner for which they were not designed, or (iv) improper storage and handling of the Products.

Current product instructions are available at <u>www.thermoscientific.com/pierce</u>. For a faxed copy, call 800-874-3723 or contact your local distributor. © 2012 Thermo Fisher Scientific Inc. All rights reserved. Unless otherwise indicated, all trademarks are property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.