



00-2213 500 ml  
PNPP (1X)-Substrate  
for Alk. Phosphatase  
Lot No.

**PNPP (p-nitrophenyl phosphate)**

**SUBSTRATE FOR ALKALINE PHOSPHATASE**

**FORM:**

Liquid (500 ml). Ready to use.

**RECOMMENDED USAGE:**

To use as a soluble substrate for alkaline phosphatase in enzyme immunoassay. Dispense the amount required for the experiment. Incubate the substrate with enzyme conjugate for 30 minutes. Read optical density at 405 nm.

**STORAGE:**

Store at 2-8°C.

**REMARKS:**

For research use only. Not for human or drug use. No preservative added.

[www.invitrogen.com](http://www.invitrogen.com)

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288

E-mail: [techsupport@invitrogen.com](mailto:techsupport@invitrogen.com)

PI002213

(Rev 10/08) DCC-08-1089

**Important Licensing Information** - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, [www.invitrogen.com](http://www.invitrogen.com)). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.



00-2213 500 ml  
PNPP (1X)-Substrate  
for Alk. Phosphatase  
Lot No.

**PNPP (p-nitrophenyl phosphate)**

**SUBSTRATE FOR ALKALINE PHOSPHATASE**

**FORM:**

Liquid (500 ml). Ready to use.

**RECOMMENDED USAGE:**

To use as a soluble substrate for alkaline phosphatase in enzyme immunoassay. Dispense the amount required for the experiment. Incubate the substrate with enzyme conjugate for 30 minutes. Read optical density at 405 nm.

**STORAGE:**

Store at 2-8°C.

**REMARKS:**

For research use only. Not for human or drug use. No preservative added.

[www.invitrogen.com](http://www.invitrogen.com)

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288

E-mail: [techsupport@invitrogen.com](mailto:techsupport@invitrogen.com)

PI002213

(Rev 10/08) DCC-08-1089

**Important Licensing Information** - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, [www.invitrogen.com](http://www.invitrogen.com)). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.