



G00-2201S 50 mg
PNPP - Substrate
for Alkaline Phosphatase
Lot No.

PNPP (p-nitrophenyl phosphate) Sample

SUBSTRATE FOR ALKALINE PHOSPHATASE

FORM: Crystalline powder (50 mg). Dissolve contents into 0.5 ml of deionized water. Mix well and keep it frozen as a stock solution. Protect from light until ready to use.

RECOMMENDED WORKING DILUTION: To use as a soluble substrate for alkaline phosphatase in an enzyme immunoassay, dilute the stock PNPP solution 1:100 with 0.75 M 2-amino-2-methyl-propanediol, pH 10.3. Discard the unused portion of the diluted PNPP. Read optical density at 405 nm using a reference filter of 490nm.

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

www.invitrogen.com

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

E-mail: techsupport@invitrogen.com

PIG002201S

(Rev 04/09) DCC-09-0417

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). 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.



G00-2201S 50 mg
PNPP - Substrate
for Alkaline Phosphatase
Lot No.

PNPP (p-nitrophenyl phosphate) Sample

SUBSTRATE FOR ALKALINE PHOSPHATASE

FORM: Crystalline powder (50 mg). Dissolve contents into 0.5 ml of deionized water. Mix well and keep it frozen as a stock solution. Protect from light until ready to use.

RECOMMENDED WORKING DILUTION: To use as a soluble substrate for alkaline phosphatase in an enzyme immunoassay, dilute the stock PNPP solution 1:100 with 0.75 M 2-amino-2-methyl-propanediol, pH 10.3. Discard the unused portion of the diluted PNPP. Read optical density at 405 nm using a reference filter of 490nm.

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

www.invitrogen.com

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

E-mail: techsupport@invitrogen.com

PIG002201S

(Rev 04/09) DCC-09-0417

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). 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.