

Protocol for Preparation of BCIP-T/NBT Substrate Solution

This protocol is for the Protocol for Preparation of BCIP-T/NBT Substrate Solution

Add the following components to 10 ml of the Alkaline Phosphatase (AP) buffer (100 mM Tris-HCl (pH 9.5), 100 mM NaCl and 10 mM MgCl₂):

BCIP-T (50 mg/ml in dimethylformamide)	33 µl
NBT (75 mg/ml in 70%dimethylformamide)	44 µl

Note

- Avoid exposure to light.
- Prepare fresh developing solution and use within an hour.
- BCIP-T and NBT solutions in dimethylformamide do not freeze. They are stable for approximately 2 years when stored at -20°C in the dark.

thermoscientific.com

© 2012 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

North America
Technical Services:
techservice.genomics@thermofisher.com
Customer Services:
customerservice.genomics@
thermofisher.com
Tel 800 235 9880
Fax 800 292 6088

Europe and Asia
Technical Services:
techservice.emea.genomics@
thermofisher.com
Customer Services:
customerservice.emea.genomics@
thermofisher.com

Thermo
S C I E N T I F I C

Part of Thermo Fisher Scientific