

PRODUCT INFORMATION

**10X Thermo Scientific Tango
Buffer** (with BSA)

_ _ _ _

Lot: _ _

Store at -20°C

BSA included

Description

Thermo Scientific Tango buffer, being a part of Five Buffer System, is specially designed for use in double digestions. For more information see www.thermoscientific.com/doubledigest.

1X Buffer Composition

33 mM Tris-acetate (pH 7.9 at 37°C),
10 mM magnesium acetate,
66 mM potassium acetate,
0.1 mg/ml BSA.

CERTIFICATE OF ANALYSIS

Functional test

Functionally tested in DNA digestion with 15 randomly picked conventional restriction endonucleases. Digestion conditions are as described in activity assays of each enzyme. Reaction products analyzed by agarose electrophoresis and ethidium bromide staining. No change in digestion pattern is observed compared to previous lot.

Endo- and Exonucleases Assay

Single stranded and double stranded radiolabeled oligonucleotides are incubated with the 1X buffer at 37°C for 16 hours, separated on polyacrylamide gel and analyzed by phosphoimaging. No signs of radiolabeled oligo degradation is considered as the absence of endo- and exonucleases activities.

Nicking activities Assay

Supercoiled pUC19 DNA is incubated with the 1X buffer at 37°C for 17 hours and analyzed by agarose gel electrophoresis and ethidium bromide staining. There is no linearised plasmid form and relaxation of supercoiled plasmid form compared to negative control.

Quality authorized by:



Jurgita Zilinskiene

Important Note

This buffer contains BSA. In order to avoid the risk of bacterial growth, do not store buffer containing BSA at 4°C for periods longer than one week. If buffer is to be stored at 4°C it should be aliquoted into small portions and the main stock should be kept at -20°C.

PRODUCT USE LIMITATION

This product is developed, designed and sold exclusively *for research purposes and in vitro use only*. The product was not tested for use in diagnostics or for drug development, nor is it suitable for administration to humans or animals.

Please refer to www.thermoscientific.com/onebio for Material Safety Data Sheet of the product.

© 2012 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.