

TNBSA

(2,4,6-Trinitrobenzene Sulfonic Acid)

TS-28997

0864w

Product Description

Number	Description
TS-28997	TNBSA (2,4,6-Trinitrobenzene Sulfonic Acid); 5% w/v in methanol, 100 ml

Introduction

2,4,6-Trinitrobenzene Sulfonic Acid (TNBSA or TNBS) is a rapid and sensitive assay reagent for the determination of free amino groups.¹ Primary amines, upon reaction with TNBSA, form a highly chromogenic derivative, which can be measured at 335 nm (see figure). Qualitative measurements of amines, sulfhydryls or hydrazides,³ and quantitative measurements of Σ -amino groups of L-lysine⁴ have also been obtained using TNBSA.

Example Protocol

The following protocol is adapted from a procedure described by Hermanson.⁴

Materials Required

Reaction Buffer: 0.1 M sodium bicarbonate, pH 8.5

TNBSA: 0.01% (w/v) solution of TNBSA. Prepare using reaction buffer as a diluent. Prepare fresh for each reaction.

10% solution of SDS in distilled water

1 N HCl

Method

1. Dissolve proteins to be assayed directly in reaction buffer at a concentration of 20-200 $\mu\text{g/ml}$. Alternatively, for proteins already in solution, the buffer can be changed by dialysis. Small molecules such as amino acids should be dissolved in reaction buffer at a concentration of 2-20 $\mu\text{g/ml}$.
Caution: Tris, glycine or other buffers containing free amines should be avoided because the free amines will react with TNBSA.
2. Add 0.25 ml of the 0.01% (w/v) solution of TNBSA to 0.5 ml of each sample solution. Mix well.
3. Incubate at 37°C for two hours.
4. Add 0.25 ml of 10% SDS and 0.125 ml of 1 N HCl to each sample.
5. Measure the absorbance of the solutions at 335 nm.

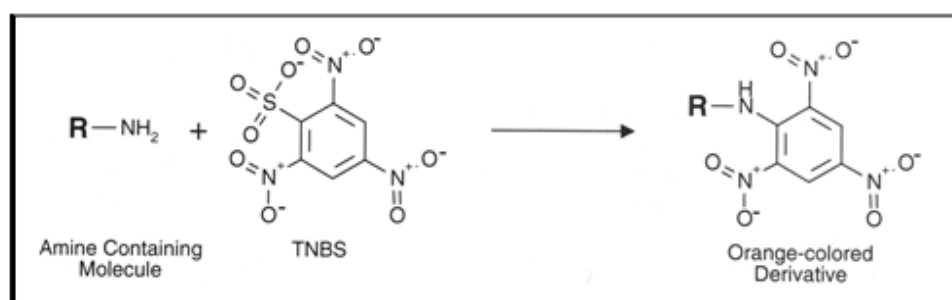
Quantitative determination of the number of amines contained within a sample can be accomplished through comparison to a standard curve generated by the use of an amine-containing compound (*e.g.*, amino acid) dissolved in a series of known concentrations. The standards should be dissolved or dialyzed into the reaction buffer and must be assayed under reaction conditions identical to those utilized for the samples.

References

1. Habeeb, A.F.S.A. (1966). Determination of free amino groups in protein by trinitrobenzene sulfonic acid. *Anal. Biochem.* **14**, 328.
2. Inman, J.K., Dintzis, H.M. (1969). *Biochem.* **8**, 4074.
3. Sashidhar, R.B., Capoor, A.K., Ramana, D. (1994). Quantitation of Σ -amino group using amino acids as reference standards by trinitrobenzene sulfonic acid. *J. Immunol. Meth.* **167**, 121-127.
4. Hermanson, G. (1996). *Bioconjugate Techniques*, p.112-113. Academic Press, San Diego, California.
This book is available from Pierce as Prod. No. 20002.
5. Endo, N., Umemoto, N., Kato, Y., Takeda, Y. and Hara, H. (1987). A novel covalent modification of antibodies at their amino groups with retention of antigen-binding activity. *J. Immunol.* **104**, 253-258.
6. Sashidhar, R.B., Capoor, A.K. and Ramana, D. (1994). Quantitation of Σ -amino group using amino acids as reference standards by trinitrobenzene sulfonic acid. *J. Immunol.* **167**, 121-127.
7. Morcol, T., Subramanian, A. and Velander, W. (1997). Dot-blot analysis of the degree of covalent modification of proteins and antibodies at amino groups. *J. Immunol.* **203**, 45-53.
8. Bubnis, W. and Ofner, C. (1992). The determination of epsilon-amino groups in soluble and poorly soluble proteinaceous materials by a spectrophotometric method using trinitrobenzenesulfonic acid. *Anal Biochem.* **297**, 129-33.

Current versions of product instructions are available at www.thermo.com/columns.

© 2008 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.



Reaction of TNBSA with a primary amine-containing molecule to produce a chromogenic derivative.