

Bovine Serum Albumin

Cat. No. 15561-020

Conc.: 50 mg/ml

Size: 150 mg

Store at -20°C (not frost-free)

Description

Bovine serum albumin (BSA) is acetylated to inactivate contaminating nucleases and proteases (1). It is suitable for use as an enzyme stabilizer during purification or for dilution of restriction endonucleases and nucleic acid modifying enzymes. BSA is also commonly used in DNA and protein labeling experiments as a blocking agent to minimize background.

Quality Control

The Certificate of Analysis provides detailed quality control information for this product. Certificates of Analysis are available on our website. Go to www.invitrogen.com/support and search for the Certificate of Analysis by product lot number, which is printed on the box.

References

1. Gonzales, N., Wiggs, J. Chamberlin, M.J. (1977), *Arch. Biochem. Biophys.* 182, 404.

©2001–2008 Invitrogen Corporation. All rights reserved.

For research use only. Not intended for any animal or human therapeutic or diagnostic use.