Glowing products for science

## Mix-n-Stain™ Total Protein Prestain Kits

Mix-n-Stain™ Total Protein Prestain Kits have been discontinued, and replaced by our VersaBlot™ Total Protein Normalization Kits.



## **Product attributes**

Conjugation CF®680T, CF®770T

Call us: 800-304-5357 Email: btinfo@biotium.com

## **Product Description**

Mix-n-Stain™ Total Protein Prestain Kits have been discontinued, and replaced by our <u>VersaBlot™ Total Protein Normalization Kits</u>. The VersaBlot™ kits include all components of the original prestain kits, plus a dye reversal agent that allows two-color NIR western detection to be performed after normalization.

This datasheet was generated on November 17, 2025 at 11:03:26 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/mix-n-stain-total-protein-prestain-kits/">https://biotium.com/product/mix-n-stain-total-protein-prestain-kits/</a>