



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

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
-------------	------------------

Product Description

Mix-n-Stain™ Total Protein Prestain Kits have been discontinued, and replaced by our [VersaBlot™ Total Protein Normalization Kits](#). 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 January 22, 2026 at 07:11:18 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/mix-n-stain-total-protein-prestain-kits/>