Glowing products for science

## Tween® 20

A polysorbate surfactant commonly used as a blocking and washing agent in Western blotting and other immunostaining assays.



## **Product attributes**

Storage Conditions Store at room temperature

CAS number 9005-64-5

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

## **Product Description**

Tween® 20 (polysorbate 20) is a non-ionic surfactant commonly used as a blocking and washing agent in western blotting and other immuno-staining assays. Tween® 20 is typically included in blocking buffers and washing buffers at a concentration of 0.05-0.1% to reduce non-specific antibody binding.

- Used in western blot blocking and washing buffers
- Used in ELISA washing buffers
- Non-ionic surfactant
- 50 mL size
- Store at room temperature

TWEEN is a registered trademark of Croda International LLC.

This datasheet was generated on November 3, 2025 at 11:31:03 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/tween-20/">https://biotium.com/product/tween-20/</a>