



Glowing products for science

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

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

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 July 31, 2026 at 03:01:05 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/tween-20/>