



Q-BRIGHT® EXTREME CHEMILUMINESCENT WESTERN BLOT KIT

SKU: QBD-60025



DESCRIPTION

Q-Bright® Extreme chemiluminescent western blot kit is our most sensitive luminol-based substrate for horseradish peroxidase (HRP) . Q-Bright® Extreme provides much greater sensitivity compared to Q-Bright® Endure and is comparable to many of the available commercial reagents providing mid- to low-femtogram levels of detection in western blot. The Western Blot Kit includes western blot wash buffer, Q-Bright® Western Blot Blocking Buffer, Anti-Mouse and Anti-Rabbit HRP conjugated secondary antibodies, and Q-Bright® Extreme Substrate and Stable Peroxide.

SPECIFICATIONS

Format	Solution
Unit Size	1 Kit
Storage Instructions	Store at 4°C. Product may ship at ambient temperature.
Origin	United States of America

For research use only. Not intended for therapeutic or diagnostic use in animals or humans.



DOCUMENTS

- [Safety Data Sheet](#)
- [Datasheet](#)

GALLERY IMAGES



For research use only. Not intended for therapeutic or diagnostic use in animals or humans.