

DAB Substrate Kit

DAB reacts with HRP in the presence of peroxide to yield an insoluble brown-colored product at locations where peroxidase-conjugated antibodies are bound to samples.



Product attributes

Call us: 800-304-5357

Storage Conditions

Store at 2 to 8 °C

Product Description

DAB (3,3'-diaminobenzidine) reacts with HRP in the presence of peroxide to yield an insoluble brown-colored product at locations where peroxidase-conjugated antibodies are bound to samples.

- Excellent substrate for immunohistochemical staining that requires the use of traditional counterstains and mounting media
- The brown precipitate is insoluble in alcohol and other organic solvents
- Kit includes 20X DAB substrate solution and 20X stabilized peroxide buffer, 15 mL each

This datasheet was generated on November 22, 2025 at 07:58:10 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/dab-substrate-kit/