



Glowing products for science

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

## 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

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 June 18, 2026 at 10:31:43 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/dab-substrate-kit/>