

EB Succinimidyl Ester

The amine reactive succinimidyl ester of EB (ethidium bromide) can be conjugated to peptides, proteins, drugs, polymeric materials and biomolecules with primary amine groups. The conjugates are expected to be essentially nonfluorescent until they are able to complex with nucleic acids, resulting in red fluorescence.



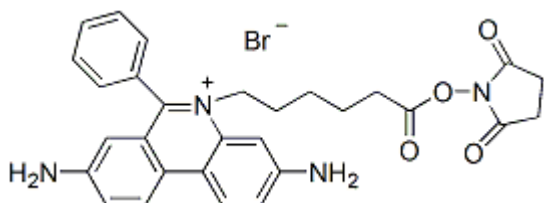
Product attributes

| Excitation/Emission | 279, 522/593 nm (with DNA) |
|---------------------|----------------------------|
|---------------------|----------------------------|

Product Description

The amine reactive succinimidyl ester of EB (ethidium bromide) can be conjugated to peptides, proteins, drugs, polymeric materials and biomolecules with primary amine groups. The conjugates are expected to be essentially nonfluorescent until they are able to complex with nucleic acids, resulting in red fluorescence. The conjugates might be useful for studies of nucleic acid binding to various biomolecules, such as DNA-binding proteins. It is also possible that conjugates of other biomolecules might be capable of monitoring their transport into the nucleus. EB dye conjugates of solid or semisolid matrices, such as microspheres, magnetic particles or various resins, might be useful for the detection or affinity isolation of nucleic acids.

- $\lambda_{Ex}/\lambda_{Em}$ (with DNA) = 522/593 nm with DNA (EB also has a large absorption peak at 279 nm)
- Red solid soluble in DMSO or DMF
- Store at -20 °C, desiccated and protect from light
- MW: 498



See our other [reactive DNA/RNA binding dyes](#).

This datasheet was generated on January 9, 2026 at 07:08:24 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/eb-succinimidyl-ester/>