

ab286890 – Safe Image Green DNA Stain

For the safe detection of dsDNA, ssDNA and RNA in agarose and polyacrylamide gels.

For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit:

<http://www.abcam.com/ab286890>

Storage and Stability

Upon arrival, the stain should be stored at 4°C. The stain is stable for 2 years from the date of shipping.

Materials Supplied

Item	Quantity	Storage Condition
Safe Image Green DNA Stain	1 mL	4°C

Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully utilize this assay:

- UV and LED transilluminators
- Pipettes
- Agarose.

Protocol

- With Safe Image Green DNA Stain, you do not need to add any dyes to the gel matrix or running buffers. Safe Image Green DNA stain is provided in a form of 6X sample loading dyes, and it alone should be added to your samples.
 - The Safe Image dyes eliminate contamination risks associated with Ethidium Bromide.
 - After the electrophoresis, view and document your results as you would do with EtBr staining protocols
1. Prepare a 100 ml agarose or polyacrylamide solution.
 2. Mix gently without introducing any air bubbles.
 3. For agarose gel, let the solution cool down to 60 - 70°C and cast the gel. For polyacrylamide gel, add APS and TEMED and cast the gel according to regular polyacrylamide gel casting protocol.
 4. Mix samples and DNA marker with Safe Image dye at a 1:5 (dye:sample) dilution rate
 5. Following electrophoresis, view the gel under UV light.

Δ Note: Safe Image Green DNA Stain is also visible under blue LED light.

Technical Support

Copyright © 2021 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

www.abcam.com/contactus

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)