## Activated Charcoal Decontamination Bags

Activated charcoal in a convenient bag for decontamination of used gel staining solutions containing ethidium bromide or other toxic dves.



## **Product attributes**

Call us: 800-304-5357

Email: btinfo@biotium.com

## **Product Description**

The Activated Charcoal Decontamination Bags were developed for convenient, no-mess clean-up of mutagenic dyes like ethidium bromide (EtBr) from used gel staining solutions.

- Remove mutagenic dyes (ie, ethidium bromide) from gel staining solutions
- Use to clean up any dyes from a solution to allow drain disposal
- No-mess, convenient bag
- Reusable bag capacity of at least 10 mg EtBr

The Activated Charcoal Decontamination Bags concentrate ethidium bromide from used gel staining solutions in a convenient bag. Concentrating the toxic dye into a small bag allows the user to dispose of it in the solid waste instead of as a large amount of liquid waste, potentially saving a lot of money in toxic waste disposal fees.

Each bag has a total capacity of at least 10 mg ethidium bromide. The bags can also be used to decontaminate other solutions containing mutagenic or toxic dyes. The Biotium gel stains GelRed® and GelGreen® have been independently certified as non-mutagenic and non-toxic, and can therefore be disposed of directly down the drain. However, since local waste procedures may vary, the Activated Charcoal Decontamination Bags can be used to remove GelRed® and GelGreen®.

The bags come with instructions for decontaminating buffers and monitoring dye removal.

GelRed and GelGreen and their uses are covered by granted and/or pending US and International patents. GelRed and GelGreen are registered trademarks of Biotium, Inc.

This datasheet was generated on November 2, 2025 at 05:53:12 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/activated-charcoal-decontamination-bags/">https://biotium.com/product/activated-charcoal-decontamination-bags/</a>