

Product Storage Statement

Product Shipping and Storage Conditions

The recommended product storage conditions can be found on the product label and in the Product Information (PI) sheet.

Most of our products are stable at room temperature for many days, but we recommend storage in a refrigerator or freezer to prevent the growth of microbes and to prolong shelf-life. Therefore, to reduce shipping costs for our customers, products with recommended storage at 4°C or -20°C may ship with an ice pack or at ambient temperature. These products may thaw without affecting product performance. When you receive the product, place it under the recommended storage conditions.

Products with recommended storage at ultra low temperature (-70°C) that also ship on dry ice should arrive frozen. If a product you received was shipped on dry ice and thawed during transit, please contact customer service.

Recommended Long-Term Storage

The storage conditions that a Biotium label recommends are deliberately conservative. Long term storage at low temperatures usually will increase the time that the product will remain within specification, unless specific instructions otherwise are provided (such as "do not refrigerate" or "do not freeze").

Temperature-Controlled Storage

Biotium labels and product information sheets may show a specific temperature for long term storage, or a temperature range that is adequate for long term storage. For all products, the implied range is as follows for the following storage conditions:

- **Room Temperature**
For Biotium products where the label indicates room temperature or RT, this implies storage in ambient conditions between 20°C and 30°C.
- **Refrigerator**
For Biotium products where the label indicates 4°C, this implies storage in a refrigerator that is between 2°C and 8°C.
- **Freezer**
For Biotium products where the label indicates -20°C, this implies storage in a freezer that is between -35°C and -5°C.
- **Ultra Low Temperature Freezer**
For Biotium products where the label indicates -70°C, this implies storage in an ultra low-temperature (ULT) freezer that is below -60°C.

Light-Sensitive Products

Fluorescent dyes should be protected from light during long-term storage. Some of our light-sensitive products are packaged in amber bottles or amber glass vials to protect the products from light. Some of our reactive dyes are packaged in transparent vials, then sealed inside a moisture-resistant foil bag, which also will protect them from light. Other fluorescent dyes in transparent vials should be stored in the dark. Fluorescent dyes generally are not sensitive to brief light exposure while they are being handled during an experiment, unless the product information sheet indicates otherwise.

Moisture-Sensitive Products

Products that are sensitive to moisture may be sealed in a moisture-resistant foil bag with a desiccant pack, which will protect them from moisture. Other products should be stored in a desiccator if indicated on the label.

Expiration Dates and Shelf Life

Assay/staining kits and PCR master mixes

The guaranteed shelf life from date of receipt for kits and master mixes is listed on the product information. Kits also may have an expiration date printed on the kit box label; this is the guaranteed shelf life calculated from the day that the product shipped from our facility.

Antibodies and other bioconjugates

The guaranteed shelf life from date of receipt for antibodies and conjugates is listed on the product information sheet.

Chemicals, dyes, and gel stains

If not otherwise specified, Biotium guarantees the stability of chemicals, dyes, and gel stains for at least one year from the date you receive the product. However, the majority of these products are highly stable, as long as they are stored as recommended.

Expired products

Products may remain functional for longer than the guaranteed shelf life. If you wish to use a Biotium product that is past its guaranteed shelf life, we recommend performing a small scale positive control experiment to confirm that the product still works for your application before processing a large number of samples or precious samples. Even if the guaranteed shelf life has expired, if any of our products fails to perform as expected, we recommend contacting technical support for troubleshooting assistance before purchasing a replacement.