



Glowing products for science

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

Water, Ultrapure Molecular Biology Grade

Suitable for use in molecular biology applications that require high quality, ultra-pure water.



Product attributes

Storage Conditions
Store at room temperature or below

Product Description

Water, Ultrapure Molecular Biology Grade is suitable for use in molecular biology applications that require high quality water.

- Molecular Biology Grade
- RNase-, DNase-, protease-, pyrogen/endotoxin-free
- No toxic chemical additives, such as DEPC
- 4L Cubitainer® with detachable spigot for easy dispensing
- Non-glug design for a continuous, uninterrupted stream
- Opening can be raised or depressed for pouring
- Not Sterile

Type I ultrapure (18 megOhm-cm) water is subjected to ultrafiltration to eliminate RNase, DNase, protease, pyrogens/endotoxins (< 0.001 EU/mL) and bacteria (< 1 cfu/mL). The product is not certified as sterile, and is therefore not recommended for cell culture. Water, Ultrapure Molecular Biology Grade is free of toxic chemical additives, such as diethyl pyrocarbonate (DEPC).

Cubitainer is a registered trademark of Hedwin Corporation.

References

1. Kartika: Jurnal Ilmiah Farmasi, 6.2, 65 (2019), [DOI: 10.26874/kjif.v6i2.153](https://doi.org/10.26874/kjif.v6i2.153)
2. Exp Ther Med.12, 4029, (2016), [DOI: 10.3892/etm.2016.3856](https://doi.org/10.3892/etm.2016.3856)
3. Life Sci, 94, 17, (2014), [DOI: 10.1016/j.lfs.2013.10.034](https://doi.org/10.1016/j.lfs.2013.10.034)

This datasheet was generated on January 8, 2026 at 04:21:13 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/water-ultrapure-molecular-biology-grade/>