



Glowing products for science

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

SuperHT PAP Pen

This item is discontinued.



Product attributes

Product Description

This item is discontinued.

References

1. Mol Genet Metab (2015) Feb;117(2):210-6 [doi:10.1016/j.ymgme.2015.11.004](https://doi.org/10.1016/j.ymgme.2015.11.004)
2. Mol Cell Proteomics (2018) Sep;17(9):1778-1787 [doi:10.1074/mcp.RA118.000680](https://doi.org/10.1074/mcp.RA118.000680)
3. Bio Protoc (2018) April 20; 8(8) [doi:10.21769/BioProtoc.2819](https://doi.org/10.21769/BioProtoc.2819)

This datasheet was generated on June 13, 2026 at 06:03:56 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/superht-pap-pen/>