

Mini Cell Scrapers

Small polyethylene cell scrapers for harvesting cells or cell lysates from 96-, 48- and 24-well plates.



Product attributes

Product Description

Mini Cell Scrapers are made of sterile plastic. Their small size is designed for harvesting cells or cell lysates from 96-, 48- and 24-well plates.

- Size: 0.5 cm (3/16") in width and 6 cm (2 3/8") in length.
- Packaged in 10 sterile packs of 20 scrapers per pack.
- Polyethylene, disposable & sterile.

References

1. Bio-protocol (2020), 10 (16) [DOI: 10.21769/BioProtoc.3716](https://doi.org/10.21769/BioProtoc.3716)
2. Front. Cell Dev. Biol. (2020) [doi: 10.3389/fcell.2020.572276](https://doi.org/10.3389/fcell.2020.572276)
3. Drug Metab Disp (2018) Feb;46(2):166-177 [doi:10.1124/dmd.117.078238](https://doi.org/10.1124/dmd.117.078238)
4. Mouse Models of Allergic Disease (2013) v1032 [10.1007/978-1-62703-496-8_7](https://doi.org/10.1007/978-1-62703-496-8_7)
5. International Journal Of Molecular Medicine (2013) 32: 1407-1414 [doi:10.3892/ijmm.2013.1514](https://doi.org/10.3892/ijmm.2013.1514)

This datasheet was generated on June 10, 2026 at 01:07:55 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/mini-cell-scrapers/>