

PRODUCT NAME TPX [®] microtubes			
Description	Cat. No.	Material	Format
1.5ml TPX [®] microtubes	M-50050	TPX [®]	50
1.5ml TPX [®] microtubes	M-50001	TPX [®]	1000
15 ml Tube TPX [®]	M-UN-15	TPX [®] (Cap: polyethylene)	50
Lot #: not available			

Product description

TPX[®] Polymethylpentene (PMP) tubes improve sonication and shearing efficiency. TPX[®] is a hard plastic which ameliorates the transmission of the ultrasound to the sample. TPX[®] tubes have been validated for use on the Bioruptor[®] family of sonication systems, including the Bioruptor[®] Standard and Bioruptor[®] Plus.

Characteristics

- Improves sonication and chromatin shearing efficiency
- Excellent transparency
- Heat Resistance

Important notice

Even though TPX[®] tubes are made of hard plastic, after prolonged or repeated sonication rounds (especially when using High Power setting), some cracks may appear in the tubes. For this reason, for sonication times exceeding 10 minutes (or for multiple 10 minutes sonication rounds), samples should be transferred to a new TPX[®] tube between sonication rounds.

It is also recommended to transfer your sample after sonication into a new tube before proceeding to the next assay steps (e.g. phenol-chloroform extraction).