

Please note: There is no manual for BK135, it is composed of three kits which have their own manuals. Check BK124-S, BK127-S and BK128-S manuals on [www.cytoskeleton.com](http://www.cytoskeleton.com) .