

**Protocol**

**Confirm Cell Lysis with NucleoBond Xtra Plasmid Purification Kits**



**Resuspend cell pellet in resuspension buffer**



**Add lysis buffer and invert 5X to mix completely**

**Cell suspension turns blue**



**Add neutralization buffer and invert to mix**

**Cell suspension starts to clear, but traces of blue remain**



**Continue mixing until the solution is homogeneous and lysis is complete**

**Cell suspension is completely colorless**

**You have now removed the SDS, renatured your plasmid DNA, and precipitated contaminating proteins and genomic DNA. Your plasmid DNA is ready to filter and use!**

Products

Cat. #	Product	Package Size	Price
740410.10	NucleoBond Xtra Midi	10 Preps	\$94.00
740410.50		50 Preps	\$442.00
740410.100		100 Preps	\$823.00
740412.10	NucleoBond Xtra Midi Plus	10 Preps	\$114.00
740412.50		50 Preps	\$523.00
740414.10	NucleoBond Xtra Maxi	10 Preps	\$217.00
740414.50		50 Preps	\$1,021.00
740414.100		100 Preps	\$1,903.00
740416.10	NucleoBond Xtra Maxi Plus	10 Preps	\$238.00
740416.50		50 Preps	\$1,141.00
740436.10	NucleoBond Xtra BAC	10 Preps	\$255.00
740436.25		25 Preps	\$573.00

Prices are subject to change without notice. Please check [www.clontech.com](http://www.clontech.com) for current pricing.

**Notice to Purchaser**

MACHEREY-NAGEL products are available to our U.S. and Canadian customers only.

Your use of these products and technologies is subject to compliance with any applicable licensing requirements described on the product's web page at <http://www.clontech.com>. It is your responsibility to review, understand and adhere to any restrictions imposed by such statements.