

4D-Nucleofector™ System - Current Software Status (01/2015)

1 4D-Nucleofector™ Standard Software

Package:	S-4D-02.16
LONZA Programs	
Version:	V02.16
Date:	20130227
Core Version:	V02.16
Serial:	***B****
X Unit Version:	V02.16
Serial:	***X****
Y Unit Version:	V02.16
Serial:	***Y****

2 4D-Nucleofector™ LogWare

Package:	S-4D-02.16
LONZA Programs	
Version:	V03.00
Date:	20130227
Core Version:	V03.00
Serial:	***B****
X Unit Version:	V03.00
Serial:	***X****
Y Unit Version:	V03.00
Serial:	***Y****