



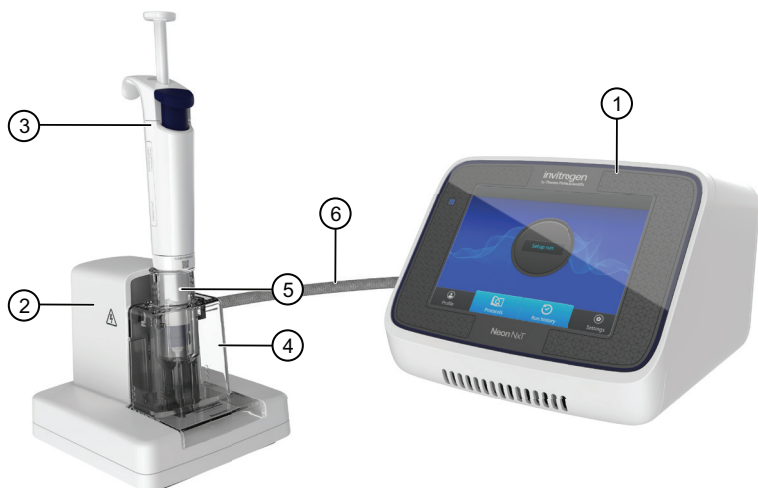
## Contents

Product	Cat. No.
Neon™ NxT Electroporation System	NEON1S
Neon™ NxT Electroporation System Starter Kit	NEON1SK



## Product description

The Neon™ NxT Electroporation System is a benchtop electroporation device that employs an electroporation technology which uses a pipette tip as an electroporation chamber to efficiently transfect mammalian cells including primary and immortalized hematopoietic cells, stem cells, and primary cells. The system efficiently delivers nucleic acids, proteins, and siRNA into all mammalian cell types including primary and stem cells with a high cell survival rate. The transfection is performed using as few as  $1 \times 10^4$  or as many as  $1 \times 10^7$  cells per reaction in a volume of 10  $\mu$ L or 100  $\mu$ L for a variety of cell culture formats (60 mm, 6-well, 48-well, 24-well, and 96-well).



### Neon™ NxT Electroporation System

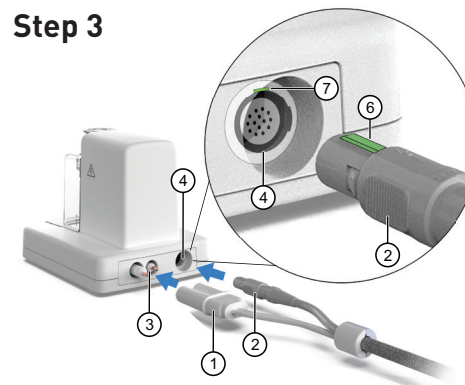
- |                              |                                    |
|------------------------------|------------------------------------|
| 1. Neon™ NxT device          | 4. Tube chamber for Neon™ NxT Tube |
| 2. Neon™ NxT Pipette Station | 5. Neon™ NxT Tube                  |
| 3. Neon™ NxT Pipette         | 6. Connector cable                 |

For Research Use Only. Not for use in diagnostic procedures.

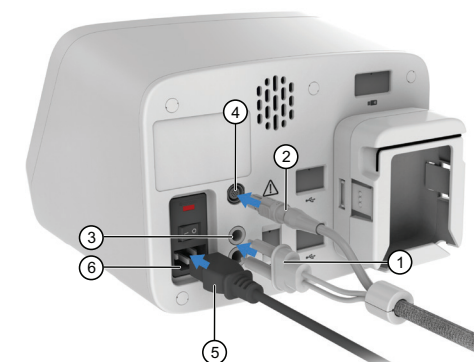
## Instrument setup

- Place the Neon™ NxT device on a level surface. Keep the area around the unit clear for proper ventilation.
- Place the Neon™ NxT Pipette Station near the Neon™ NxT device.
- Connect the high voltage ① and low voltage ② interface connectors on the cable to the high voltage ③ and low voltage ④ interface ports at the rear of the Neon™ NxT Pipette Station. **Note:** Ensure the notch ⑥ on the low voltage connector ② is aligned with the corresponding slot ⑦ on the low voltage port ④.
- Connect the high voltage ① and low voltage ② interface connectors on the cable to the high voltage ③ and low voltage ④ interface ports at the rear of the Neon™ NxT device. **Note:** Ensure the notch on the low voltage connector is properly aligned with the low voltage port.
- Ensure the AC power switch is in the **Off** position.
- Attach the power cord ⑤ to the AC inlet ⑥ at the rear of the Neon™ NxT device, then plug the cord into an electrical outlet.

### Step 3








### Step 4–6





## Online resources

- Scan the QR code for videos on how to use the Neon™ NxT pipette, set up the system, and other FAQs.
- Visit [thermofisher.com/neonnxT](https://thermofisher.com/neonnxT) for manuals, safety, and additional product information.
- For support, visit [thermofisher.com/support](https://thermofisher.com/support).



Symbol	Action
<b>Calculate electroporation volumes</b>	
 <b>Cell count calculator</b>	a. Select <b>Cell count calculator</b> b. Enter the experimental parameters c. Select <b>Calculate</b>
<b>Create a protocol</b>	
 <b>Quick run</b>	a. Select <b>Set up run &gt; Quick run</b> b. Enter the electroporation parameters c. Select <b>Save protocol</b>
<b>Create a plate map</b>	
 <b>Create plate</b>	a. Select <b>Set up run &gt; Create plate</b> b. Select plate type (6, 12, 24, or 96-well) c. Select run order (top to bottom/by order of increasing voltage) d. Select wells and assign protocols e. Select <b>Save plate</b>
<b>Edit a plate map</b>	
 <b>Open plate</b>	a. Select <b>Set up run &gt; Open plate</b> b. Select the protocol to modify c. Select wells and assign protocols d. Select <b>Save plate</b>
<b>Open a protocol or plate map</b>	
 <b>Library</b>	a. Select <b>Library &gt; Protocol Library</b> or <b>Plate Library</b> b. Select a protocol or plate map

Symbol	Function		
<b>Open Sign in screen</b>			
 <b>Sign in</b>	Select <b>Sign in</b> to log in to the instrument Provides access to the following items: <ul style="list-style-type: none"> <li>▪ Create user profile</li> <li>▪ Change PIN</li> <li>▪ Manage user profiles (Administrator only)</li> </ul>		
<b>Open Settings screen</b>			
 <b>Settings</b>	Select <b>Settings</b> Provides access to the following items: <table border="0" style="width: 100%;"> <tr> <td style="vertical-align: top;"> <ul style="list-style-type: none"> <li>▪ About instrument               <ul style="list-style-type: none"> <li>▪ EULA</li> <li>▪ Check updates</li> </ul> </li> <li>▪ Instrument settings               <ul style="list-style-type: none"> <li>▪ Instrument name</li> <li>▪ Date &amp; time</li> <li>▪ Sleep mode</li> <li>▪ Brightness</li> <li>▪ Network configuration</li> <li>▪ Cloud region</li> <li>▪ Auto sign out</li> </ul> </li> </ul> </td> <td style="vertical-align: top;"> <ul style="list-style-type: none"> <li>▪ Maintenance &amp; services               <ul style="list-style-type: none"> <li>▪ Software update</li> <li>▪ Self verification test</li> <li>▪ Export instrument log</li> <li>▪ Restore factory settings</li> </ul> </li> <li>▪ Run history               <ul style="list-style-type: none"> <li>▪ Export run report</li> <li>▪ Delete run report</li> </ul> </li> </ul> </td> </tr> </table>	<ul style="list-style-type: none"> <li>▪ About instrument               <ul style="list-style-type: none"> <li>▪ EULA</li> <li>▪ Check updates</li> </ul> </li> <li>▪ Instrument settings               <ul style="list-style-type: none"> <li>▪ Instrument name</li> <li>▪ Date &amp; time</li> <li>▪ Sleep mode</li> <li>▪ Brightness</li> <li>▪ Network configuration</li> <li>▪ Cloud region</li> <li>▪ Auto sign out</li> </ul> </li> </ul>	<ul style="list-style-type: none"> <li>▪ Maintenance &amp; services               <ul style="list-style-type: none"> <li>▪ Software update</li> <li>▪ Self verification test</li> <li>▪ Export instrument log</li> <li>▪ Restore factory settings</li> </ul> </li> <li>▪ Run history               <ul style="list-style-type: none"> <li>▪ Export run report</li> <li>▪ Delete run report</li> </ul> </li> </ul>
<ul style="list-style-type: none"> <li>▪ About instrument               <ul style="list-style-type: none"> <li>▪ EULA</li> <li>▪ Check updates</li> </ul> </li> <li>▪ Instrument settings               <ul style="list-style-type: none"> <li>▪ Instrument name</li> <li>▪ Date &amp; time</li> <li>▪ Sleep mode</li> <li>▪ Brightness</li> <li>▪ Network configuration</li> <li>▪ Cloud region</li> <li>▪ Auto sign out</li> </ul> </li> </ul>	<ul style="list-style-type: none"> <li>▪ Maintenance &amp; services               <ul style="list-style-type: none"> <li>▪ Software update</li> <li>▪ Self verification test</li> <li>▪ Export instrument log</li> <li>▪ Restore factory settings</li> </ul> </li> <li>▪ Run history               <ul style="list-style-type: none"> <li>▪ Export run report</li> <li>▪ Delete run report</li> </ul> </li> </ul>		

**Limited product warranty:** Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at [www.thermofisher.com/us/en/home/global/terms-and-conditions.html](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html).

**Disclaimer:** TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

**Manufacturer:** Life Technologies Holdings Pte Ltd Block 33 | Marsiling Industrial Estate Road 3 | #07-06, Singapore 739256

©2023 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.