

# GeneMapper™ Software v6.0 Oracle™ security patch for TNS listener vulnerability

Pub. No. MAN0024762 Rev. A.0

## Overview

This user bulletin explains how to download and install a security patch for the Oracle™ software that addresses TNS listener vulnerability.

## Download and run the RunPatch.bat utility

1. Close the GeneMapper™ Software.
2. Log in to the GeneMapper™ Software computer with administrative privileges.
3. Download the update from this link <https://downloads.thermofisher.com/GeneMapper/13329696.zip>.
4. Unzip the downloaded file: right-click on **13329696.zip**, then select **Extract All**.
5. Copy the extracted **13329696** folder to the C:\ drive.

---

**IMPORTANT!** Save the file in the root directory, *not* the desktop. The final folder path must be **C:\13329696**.

---

6. In Windows™ Explorer, open the **C:\13329696** folder.
7. Right-click **runPatch.bat**, then select **Run as administrator**.  
The utility takes ~10 minutes to run. When the utility is finished, a message is displayed: **opatch script stated...OPatch succeeded**.

---

**IMPORTANT!** If the utility fails at any time, restart the computer before you run the utility again.

---



Life Technologies Corporation | 200 Oyster Point Blvd | South San Francisco, CA 94080 | USA  
For descriptions of symbols on product labels or product documents, go to [thermofisher.com/symbols-definition](https://www.thermofisher.com/symbols-definition).

The information in this guide is subject to change without notice.

**DISCLAIMER:** TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. 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.

**Revision history:** Pub. No. MAN0024762

Revision	Date	Description
A.0	27 January 2021	User bulletin for GeneMapper v6 Oracle patch.

**Important Licensing Information:** These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

**NOTICE TO PURCHASER: DISCLAIMER OF LICENSE:** Purchase of this software product alone does not imply any license under any process, instrument or other apparatus, system, composition, reagent or kit rights under patent claims owned or otherwise controlled by Thermo Fisher Scientific, either expressly, or by estoppel.

©2021 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.