

Customer application note program

We are proud of our large collection of Application and Technical Notes, but know that our customers use our equipment, columns, software, and sample preparation products for applications we never imagined. Now our customers have the opportunity to publish their novel applications on our website as Customer Application Notes. Similar to our Application and Technical Notes, a Customer Application Note will cite the authors by name, have an assigned number (e.g., Customer Application Note 105), and have a month and year of publication. Once finalized, other customers will be able to download it from Thermoscientific.com/dionex as a PDF. If you would like to participate in this program, your submission will circulate through the following steps:

1. Customer e-mails a well-written application as a Microsoft® Word document to jeff.rohrer@thermofisher.com. Figures should be provided in either Microsoft PowerPoint (.pptx) or Adobe® Illustrator (.eps) format. If the figure contains a chromatogram, please include a Thermo Scientific Dionex Chromeleon (.cmb) file containing the chromatogram. It may be possible for us to work with other file formats, but please contact us before submitting them. The Customer Application template is somewhat flexible, but we suggest including: Title, Authors, Introduction, Experimental (which includes Equipment, Materials, and Methods), Results, and Discussion (which includes a Conclusion or a summary paragraph).
2. We will review the document for clarity, Thermo Scientific product accuracy, trademarks, and novelty relative to other Thermo Scientific Application, Technical, and Customer Application Notes. If we believe the document will make a good Customer Application Note, the document will be returned with suggested edits. We may also write a paragraph to add to the prospective Customer Application Note. This paragraph will be labeled, "Thermo Fisher Scientific Comments" and may include a variety of information, including a comparison between the Customer Application Note and existing Thermo Scientific Application, Technical, or Customer Application Notes and/or alternative Thermo Scientific equipment and consumables which can be used to update or improve the application.
3. If you wish to proceed with your submission, please implement the suggested edits, review and either edit or approve the Thermo Fisher Scientific Comments section, and return the edited and approved document along with a completed copyright form. Briefly, this form states that the figures have not been used in other publications. It also states that we will freely share the Customer Application Note figures with other parties (see below).
4. We will then format your Customer Application Note into our template and you will have a chance to review it again before it is published onto our website.

Customer Application Note Permission Form

1. I am/we are the sole author(s) of the Customer Application Note identified below, which has been submitted to Thermo Fisher Scientific Inc. ("Thermo Fisher") for publication:

2. I/we (or my employer, per section 4) hereby grant Thermo Fisher a perpetual, royalty-free worldwide nonexclusive license, under any copyright or other intellectual property right owned by me/us, to display, copy, modify, distribute and transmit the Customer Application Note in any form and for any reason whatsoever (including, without limitation, in advertising, sales literature and/or other materials promoting the products and services of Thermo Fisher).
3. Except as otherwise provided below, I/we have full rights and authority to grant the license set forth in Section 2, and have not transferred or assigned our copyright in the Customer Application Note or any portion thereof to any other party.
4. If the Customer Application Note was created within the scope of my/our employment, I/we hereby certify that the person signing below on behalf of my/our employer is empowered to grant the license set forth in Section 2.
5. I/we certify that the Customer Application Note does not disclose or contain proprietary information belonging to my employer or any other party, and that, to the best of my/our knowledge, the method or technique described therein does not infringe the intellectual property rights of any party.
6. I/we understand that Thermo Fisher owns the copyright to any modifications made by Thermo Fisher to the Customer Application note.
7. I/we understand that Thermo Fisher is not under any obligation to publish the Customer Application Note.

Author 1 _____ Title/Position _____

Signature _____ Date _____

Phone (____) _____ E-mail _____

Author 2 _____ Title/Position _____

Signature _____ Date _____

Phone (____) _____ E-mail _____

Company/Institution _____

Address _____

City _____ State/Province _____ Zip/Postal Code _____

Employer's Signature _____ Title/Position _____

Date _____

Please scan and e-mail this completed, signed document to Jeff Rohrer, Applications Development Director, at jeff.rohrer@thermofisher.com, or mail to him at Thermo Fisher Scientific, P.O. Box 3603, Sunnyvale, CA 94088-3603, USA.