


Axiom™ Propel XPRES Reagent Kit, 2x384HT

Catalog Numbers 952352

Pub. No. MAN0018847 Rev. A.0

 **WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from thermofisher.com/support.

Product use

The Applied Biosystems™ Axiom™ Propel XPRES Reagent Kit, 2x384HT enables genotyping of SNPs and insertions/deletions (indels) (on Axiom™ genome-wide array plates and Axiom™ myDesign™ Array Plates). The kits include the reagents that are required to:

- Produce hybridization-ready targets for Axiom™ genotyping array plates, Axiom™ myDesign™ Array Plates using the Thermo Scientific™ Multidrop™ Combi Reagent Dispenser.
- Process array plates on the Applied Biosystems™ GeneTitan™ Multi-Channel Instrument.

The Axiom™ Propel XPRES Reagent Kit, 2x384HT includes Fast Target Preparation reagents which decrease the amplification time to 4 hours. Fast GeneTitan™ Reagents offer 1+ hours faster array washing and staining. Contact Thermo Fisher Scientific support to obtain the necessary library files before conducting array processing on the GeneTitan™ Multi-Channel Instrument.

Ordering information

Cat. No.	Product name	Description
952352	Axiom™ Propel XPRES Reagent Kit, 2x384HT	Sufficient for two 384HT array format plates. Includes: <ul style="list-style-type: none"> • Axiom™ Propel XPRES Reagent Kit Module 1 for 384HT only • Axiom™ Propel Reagent Kit Module 2-1 for 96F or 384HT • Axiom™ Propel Reagent Kit Module 2-2 for 96F or 384HT • Module 3 • Axiom™ Propel Fast Wash Reagent Kit Module 4-1 for 96F or 384HT • Axiom™ Propel Reagent Kit Module 4-2 for 96F or 384HT
952392	Axiom™ Propel Target Preparation Reagents, 2x384HT	Sufficient for two 384HT array format plates. Includes: <ul style="list-style-type: none"> • Axiom™ Propel XPRES Reagent Kit Module 1 for 384HT only • Axiom™ Propel Reagent Kit Module 2-1 for 96F or 384HT • Axiom™ Propel Reagent Kit Module 2-2 for 96F or 384HT
952393	Axiom™ Propel GeneTitan™ Fast Reagents, 2x384HT	Sufficient for two 384HT array format plates. Includes: <ul style="list-style-type: none"> • Module 3 • Axiom™ Propel Fast Wash Reagent Kit Module 4-1 for 96F or 384HT • Axiom™ Propel Reagent Kit Module 4-2 for 96F or 384HT

Instructions for use

For instructions on using this reagent kit, see the *Axiom™ Propel XPRES 384HT Workflow User Guide*, Pub. No. MAN0018847.

Contents and storage

IMPORTANT! The Applied Biosystems™ Axiom™ Propel XPRES Reagent Kit, 2x384HT is for single use only. Large fill reagent kits are configured to include priming volumes for the Multidrop™ Combi cassettes and have been incorporated into the master mix formulations. Discard all excess reagents after use.

Component	Storage
Axiom™ Propel XPRES Reagent Kit Module 1 for 384HT only Axiom™ Propel 10X Denat Solution Axiom™ Propel Neutral Solution Axiom™ Propel Water Axiom™ Propel Amp Solution Axiom™ XPRES Amp Enzyme	-25°C to -15°C
Axiom™ Propel Reagent Kit Module 2-1 for 96F or 384HT Axiom™ Propel Frag Enzyme Axiom™ Propel 10X Frag Buffer Axiom™ Propel Precip Solution 2 Axiom™ Propel Hyb Buffer Axiom™ Propel Hyb Solution 1	-25°C to -15°C
Axiom™ Propel Reagent Kit Module 2-2 for 96F or 384HT Axiom™ Propel Frag Diluent Axiom™ Propel Frag Reaction Stop Axiom™ Propel Precip Solution 1 Axiom™ Propel Resuspension Buffer Axiom™ Propel Hyb Solution 2	2°C to 8°C
Module 3 Axiom™ Wash Buffer A Axiom™ Wash Buffer B Axiom™ Water	Room temperature
Axiom™ Propel Fast Wash Reagent Kit Module 4-1 for 96F or 384HT Axiom™ Propel Ligation Buffer Axiom™ Fast Ligation Enzyme Axiom™ Propel Ligation Solution 1 Axiom™ Propel Probe Mix 1 Axiom™ Propel Stain Buffer Axiom™ Propel Stabilize Solution	-25°C to -15°C
Axiom™ Propel Reagent Kit Module 4-2 for 96F or 384HT Axiom™ Propel Wash A Axiom™ Propel Probe Mix 2 Axiom™ Propel Ligation Solution 2 Axiom™ Propel Stain 1-A Axiom™ Propel Stain 2-A Axiom™ Propel Stabilize Diluent Axiom™ Water Axiom™ Propel Hold Buffer Axiom™ Propel Stain 1-B Axiom™ Propel Stain 2-B	2°C to 8°C

Component part numbers

Note: The following component part numbers are for identification purposes only and cannot be used to order components separately from the reagent kits.

Component	Part No.
Axiom™ Propel XPRES Reagent Kit Module 1 for 384HT only	952351
Axiom™ Propel 10X Denat Solution	952176
Axiom™ Propel Neutral Solution	952173
Axiom™ Propel Water	952177
Axiom™ Propel Amp Solution	952174
Axiom™ XPRES Amp Enzyme	952349
Axiom™ Propel Reagent Kit Module 2-1 for 96F or 384HT	952263
Axiom™ Propel Frag Enzyme	952181
Axiom™ Propel 10X Frag Buffer	952179
Axiom™ Propel Precip Solution 2	952178
Axiom™ Propel Hyb Buffer	952182
Axiom™ Propel Hyb Solution 1	952183
Axiom™ Propel Reagent Kit Module 2-2 for 96F or 384HT	952265
Axiom™ Propel Frag Diluent	952184
Axiom™ Propel Frag Reaction Stop	952190
Axiom™ Propel Precip Solution 1	952203
Axiom™ Propel Resuspension Buffer	952206
Axiom™ Propel Hyb Solution 2	951979
Module 3	—
Axiom™ Wash Buffer A	901446
Axiom™ Wash Buffer B	901447
Axiom™ Water	901578
Axiom™ Propel Fast Wash Reagent Kit Module 4-1 for 96F or 384HT	952369
Axiom™ Propel Ligation Buffer	952208
Axiom™ Fast Ligation Enzyme	952367
Axiom™ Propel Ligation Solution 1	952212
Axiom™ Propel Probe Mix 1	952213
Axiom™ Propel Stain Buffer	952214
Axiom™ Propel Stabilize Solution	952215
Axiom™ Propel Reagent Kit Module 4-2 for 96F or 384HT	952268
Axiom™ Propel Wash A	952218
Axiom™ Propel Probe Mix 2	952217
Axiom™ Propel Ligation Solution 2	952216
Axiom™ Propel Stain 1-A	952219
Axiom™ Propel Stain 2-A	952231
Axiom™ Propel Stabilize Diluent	952248
Axiom™ Water	952177
Axiom™ Propel Hold Buffer	952254
Axiom™ Propel Stain 1-B	952258
Axiom™ Propel Stain 2-B	952260

Related documentation

- *GeneTitan™ MC Protocol for Axiom™ Array Plate Processing Quick Reference* (Pub. No. MAN0017718)

Customer and technical support

Visit [thermofisher.com/support](https://www.thermofisher.com/support) for the latest service and support information.

- Worldwide contact telephone numbers
- Product support information
 - Product FAQs
 - Software, patches, and updates
 - Training for many applications and instruments
- Order and web support

- Product documentation
 - User guides, manuals, and protocols
 - Certificates of Analysis
 - Safety Data Sheets (SDSs; also known as MSDSs)

Note: For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

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 at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



Thermo Fisher Scientific Baltics UAB | V.A. Graiciuno 8, LT-02241 | Vilnius, Lithuania
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. MAN0018847

Revision	Date	Description
A.0	09 July 2020	New publication.

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.

©2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.