

# TrueDesign Genome Editor

Workflow guide for using the TrueTag Knockout Enrichment kit for knockout cell line generation

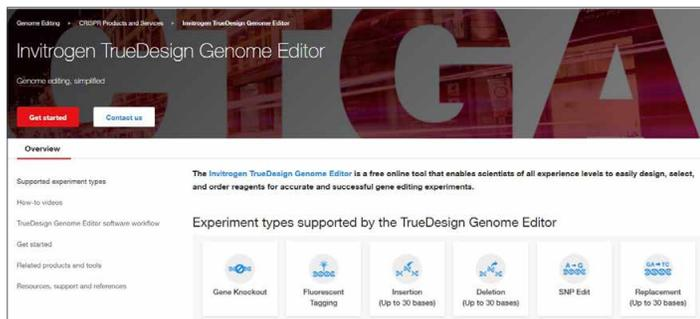
The Invitrogen™ TrueDesign™ Genome Editor is easy-to-use, free online software for designing and ordering the reagents needed for precise genome editing by homology-directed repair with RNA-guided nucleases and single-stranded DNA donors.

This workflow guide will walk through the steps for creating a gene knockout and enabling enrichment of edited cells with the Invitrogen™ TrueTag™ Knockout Enrichment Donor DNA Kit. This kit provides a template for insertion of a cassette with stop codons, GFP/RFP reporters, and blasticidin and puromycin resistance markers.

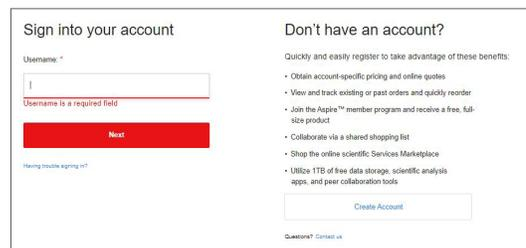
Please note that algorithm-designed CRISPR gRNAs optimized for functional knockout are available for all human and mouse genes and may be ordered directly from [thermofisher.com/truedesign](https://thermofisher.com/truedesign).

## Step 1:

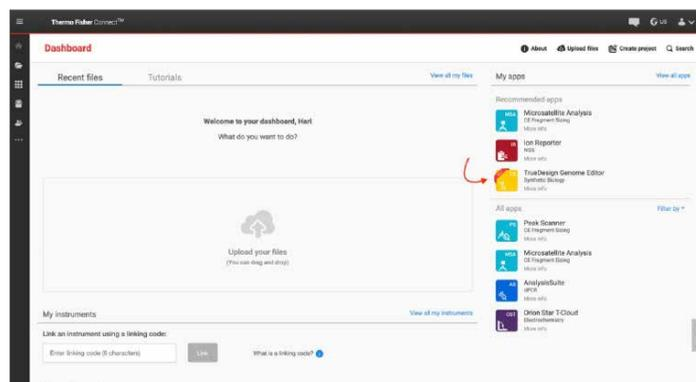
Go to [thermofisher.com/truedesign](https://thermofisher.com/truedesign). Select one of the links to launch the software.



You may be prompted to sign in. Use your existing credentials, or simply provide an email address to register as a new user.



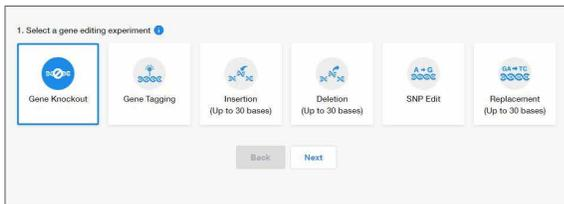
Alternatively, go directly to the **Thermo Fisher™ Connect Platform** and navigate to the TrueDesign Genome Editor.



A “Terms of Use” window may pop up. Read the content, scroll to the bottom of the screen, and click “Accept.”

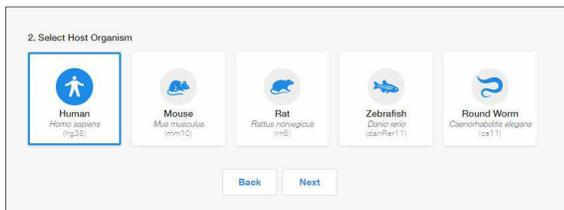
## Step 2:

In the TrueDesign software, choose **Gene Knockout** as your experiment type and click “Next.”



## Step 3:

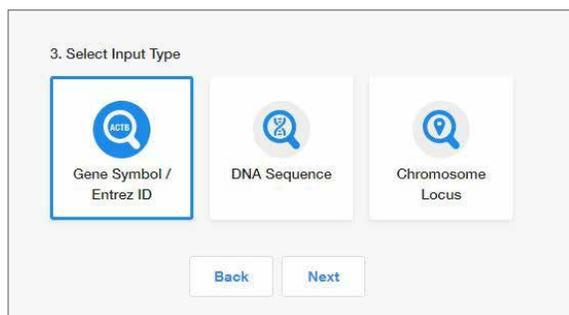
Select the host organism for your knockout experiment and click “Next.”



## Step 4:

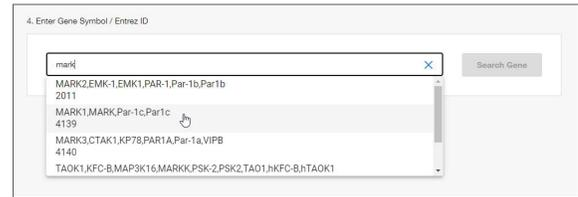
Select **Gene Symbol/Entrez ID** to identify your gene of interest. You may also navigate to a transcript of interest by entering a DNA sequence or chromosome location.

Click “Next.”



## Step 5:

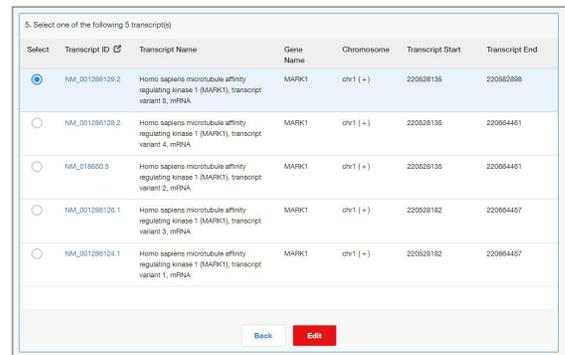
Begin typing the gene symbol or Entrez ID in the gene identifier box. A filtered drop-down list will appear. Select your gene of interest and click “Search Gene.”



## Step 6:

All of the protein-coding transcripts for your selected gene will be displayed. If there is more than one protein-coding transcript and you are unsure of which one to select, click the transcript ID hyperlink to be taken to the NCBI website, where you can better visualize the transcript maps.

After you make a selection, click “Edit.”



Select	Transcript ID	Transcript Name	Gene Name	Chromosome	Transcript Start	Transcript End
<input checked="" type="radio"/>	<a href="#">NM_001286128.2</a>	Homo sapiens microtubule affinity regulating kinase 1 (MARK1), transcript variant 5, mRNA	MARK1	chr1 (-)	220528135	220582698
<input type="radio"/>	<a href="#">NM_001286128.2</a>	Homo sapiens microtubule affinity regulating kinase 1 (MARK1), transcript variant 4, mRNA	MARK1	chr1 (+)	220528135	220664461
<input type="radio"/>	<a href="#">NM_018800.5</a>	Homo sapiens microtubule affinity regulating kinase 1 (MARK1), transcript variant 2, mRNA	MARK1	chr1 (+)	220528135	220664461
<input type="radio"/>	<a href="#">NM_001286126.1</a>	Homo sapiens microtubule affinity regulating kinase 1 (MARK1), transcript variant 3, mRNA	MARK1	chr1 (+)	220528182	220664457
<input type="radio"/>	<a href="#">NM_001286124.1</a>	Homo sapiens microtubule affinity regulating kinase 1 (MARK1), transcript variant 1, mRNA	MARK1	chr1 (+)	220528182	220664457

## Step 7:

The next screen will display the entire transcript’s topology along the top of the screen, with a zoomed-in sequence-level view below.

You will be prompted to choose **Frameshift Indels**, **Insert STOP Codon**, or **Knockout Enrichment**.

- **Frameshift indel:** The most common approach for gene knockout, this process relies on imperfect repair following a double-stranded DNA break. You will be required to select a region on the transcript to search for custom CRISPR and TALEN™ designs. Predesigned gRNAs for human and mouse are available with the **Invitrogen™ TrueGuide™ Synthetic gRNA** product line.



## Step 10 (continued)

To view the TALEN targets, click on the TALEN tab of the table, and similar information will be displayed for each TALEN pair. TAL effector nuclease (TALEN) pairs are recommended when there are no PAM sites within 10 bp of the knock-in site, or if the efficiency and specificity of the gRNAs are not optimal. Green checkmarks in the design results table will indicate the recommended technology. Learn more about TALEN technology at [thermofisher.com/tal](http://thermofisher.com/tal).

## Step 11:

To select one or more CRISPR gRNAs or TALEN pairs to add to your experiment, use the checkboxes in the table and click “Next.”

## Step 12:

The summary page will display all the reagents needed for your knockout enrichment experiment and give you the opportunity to add additional products to complete your workflow.

The screenshot displays the 'Reagents for your Gene Knockout experiment' interface. It is divided into several sections:

- CRISPR Reagents:** Includes checkboxes for CRISPR gRNA Format, Cas9 Format, Cas9 Reagent, Transfection Reagent, and TrueTag Donor Primers. Each has a dropdown menu for selection.
- TALEN Reagents:** Includes checkboxes for TALEN Format, Transfection Reagent, and TrueTag Donor Primers.
- GCD/Sequencing Primers:** Includes a checkbox for Primers and a dropdown for GCD Primer pairs.
- Reagents for Knockout Enrichment experiment:** Includes checkboxes for TrueTag Knockout Enrichment Donor DNA Kit and TrueTag Verification Primers.
- Other Products You May Need:** A table listing additional products and their quantities.

Product Name	Quantity
<input type="checkbox"/> TrueGuide sgRNA Positive Control AAVS1 (Human)	3.0 nmol * 1
<input type="checkbox"/> TrueGuide sgRNA Positive Control CDK4 (Human)	3.0 nmol * 1
<input type="checkbox"/> TrueGuide sgRNA Positive Control HPRT1 (Human)	3.0 nmol * 1
<input type="checkbox"/> TrueGuide sgRNA Negative Control	3.0 nmol * 1
<input type="checkbox"/> GeneArt Genomic Cleavage Detection Kit	1 Kit

At the bottom, there are buttons for 'Download Designs and Protocol', 'Custom Services', 'Back', 'Start a new experiment', and 'Add to Cart'.

For a knockout enrichment experiment using CRISPR technology, the gRNAs and Invitrogen™ TrueTag™ donor primers selected in the design step are automatically included. The tool will also add the TrueTag Knockout Enrichment Donor DNA Kit, which includes the required donor templates, PCR reagents, and cleanup kit to generate your transfection-ready donor DNA. Also included are Invitrogen™ TrueTag™ verification primers to do junction analysis to ensure proper insertion of the donor template.

Products may be deselected from the product summary area, or different quantities may be selected from the drop-down lists.

## Step 13:

You may use the checkboxes to add additional items such as sequencing primers or positive and negative experimental controls.

When you have completed your product selections, click “Add to Cart” for easy one-step ordering of all selected reagents.

If “Add to Cart” is not enabled in your region or you want to send the list of reagents to your purchasing agent, you can download and save a detailed report of your experiment by clicking “Download Designs.” The resulting Microsoft™ Excel™ file contains multiple tabs that include all the designs generated by the software, plus all of the gene-specific experimental details and ordering information.



Get started at [thermofisher.com/truedesign](http://thermofisher.com/truedesign)

**For Research Use Only. Not for use in diagnostic procedures.** © 2021, 2022 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Microsoft and Excel are trademarks of Microsoft Corporation. TALEN is a trademark of Collectis Bioresearch. **COL117781 0322**

**ThermoFisher**  
SCIENTIFIC