



## MMDx Kidney

# The Molecular Microscope<sup>®</sup> Diagnostic System (MMDx<sup>®</sup>) for Kidney Transplant Patients

As part of your transplant journey, your doctor may recommend a biopsy to assess the health of your kidney. A biopsy will provide information that your clinical team can use to guide your post-transplant treatment and to detect any signs of rejection.

A biopsy is typically performed as an out-patient procedure with local anesthetic. During the procedure, a very small sample is taken from your kidney.

Based on a decade of research, MMDx Kidney is a cutting-edge test for transplant biopsies. It will help your clinical team make an accurate diagnosis and smart decisions about your treatment.

### What is MMDx Kidney?

When your doctor performs a biopsy, they may use **MMDx Kidney, a test designed to assess the health of your transplant.**

The process is very straightforward. Your biopsy sample is sent to a reference lab, where the MMDx Kidney test is used to measure the expression of specific genes in your transplanted organ.

If your transplant is injured or being damaged by your immune system, specific genes are “expressed” by the kidney: information encoded in DNA is transcribed by the cell to mRNA (messenger RNA).

By detecting and reading the patterns in the mRNA, **MMDx Kidney can recognize rejection or injury** in the molecules of your kidney. This information may help your doctor to determine if you need treatment and, if necessary, quickly and accurately individualize your treatment options.

## How does this help me?

This test helps your clinical team by providing additional information regarding the health of your transplanted kidney.

MMDx Kidney can distinguish rejection or injury at the molecular level. If other diagnostic tests have proved inconclusive, MMDx Kidney may help your doctor determine whether you need additional treatment. If you do, the precise and objective information provided by **MMDx Kidney can help your doctor recommend a treatment plan that's personalized for you.**

## Does this mean I need another biopsy?

No, this can be done with your scheduled biopsy, which is either for surveillance or is requested by your doctor if your kidney is not functioning as expected.

## What will this cost me?

Like any diagnostic procedure, the cost of the test may vary based on your insurance and/or healthcare provider.

If you're interested in trying MMDx Kidney for your next biopsy, ask your doctor to inquire about the cost. The dedicated team at the MMDx reference lab will support your doctor in the prior authorization and billing appeal process.

## Is MMDx covered by my insurance provider?

A growing number of insurance companies are reimbursing the cost of MMDx Kidney, including Medicare. The lab that runs the MMDx Kidney test has a dedicated team who works closely with patients, physicians, and insurance companies to obtain coverage through accurate billing requests and appeals.

## How do I get MMDx?

Ask your physician if MMDx Kidney is appropriate for your next biopsy. A specimen kit can be ordered at [thermofisher.com/onelambda/mmdx](https://thermofisher.com/onelambda/mmdx).

## How quickly will I get my results?

Your doctor will receive your test results within 48 hours after the reference lab receives your biopsy sample.

Learn more at [go.1lambda.com/mmdxkidney](https://go.1lambda.com/mmdxkidney)  
or email us at [1lambda-mmdx@thermofisher.com](mailto:1lambda-mmdx@thermofisher.com)

MMDx and Molecular Microscope are registered trademarks of Transcriptome Sciences Inc.

MMDx Heart and MMDx Kidney are tests that were developed and validated by Kashi Clinical Laboratories, Inc. The laboratory developed tests are used for clinical purposes by the CLIA-certified laboratory performing the test. These tests have not been cleared or approved by the US FDA or CE marked in the EU as an in vitro diagnostic test.

© 2022 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.