

Electronic Certificate of Analysis Documents

Promega Corporation
2800 Woods Hollow Road, Madison, WI 53711



This project saves over a half million sheets of paper—over 1,000 reams of paper—not to mention the ink toner and electricity needed to produce these documents.

Introduction

Promega is committed to minimizing materials used and waste generated at all steps of the product lifecycle. This includes supporting documents that are shipped with kits, like a Certificate of Analysis. This document is issued by our Quality Assurance department to confirm that a specific lot of a product has met the product specifications. These documents are currently available on **Promega.com** in an easily searchable database. Historically, each product that has an issued Certificate of Analysis has also included a paper copy with the shipment.

Project Description

To support our sustainability aspirations, Promega will no longer ship paper copies of Certificates of Analysis for more than 1,000 products. This will save over 500,000 sheets of paper every year. Continuing to provide these documents electronically, rather than in hard copy, provides customers with easy-to-access product quality information and reduces the amount of paper used by everyone. Even with recycling efforts to reduce the environmental effect of paper usage, consuming less paper is a more effective way to minimize our use of limited natural resources.

Reducing our paper consumption saves trees, energy and water, and avoids pollution. Environmental benefits from going paperless are recognized at all aspects of the paper lifecycle—raw material extraction, production, transportation and disposal. Trees play a critical role in sequestering carbon, and provide oxygen to support life on our planet. Moving to electronic documents is just one way that we are conserving our natural resources.

Learn more about our sustainability efforts at:

www.promega.com/responsibility

Learn More

Supporting electronic document efforts helps us all reduce paper usage. If you are looking for a Certificate of Analysis, there are two easy ways to find them—from the **Certificate of Analysis** page, or from the product page itself.

Questions? Contact our Technical Services team at: techserv@promega.com