



Revvity Transcribe AI

Automating efficiency.

Transform your labs workflow

Tired of manually entering data from Dry Blood Spot (DBS) cards? **Revvity Transcribe AI** streamlines your process by converting handwritten text from scanned DBS card images into a digital format, speeding up processing and ensuring data accuracy—so you can focus on what matters most.

Key benefits


- **Speeds up processing time:** Digitize DBS cards faster than ever.
- **AI-powered conversion:** Transforms handwritten text into digital format with ease.
- **Minimize error:** Designed to reduce errors in data transcription.
- **Flexible** and designed to integrate with lab workflows easily.

How Revvity Transcribe AI works




Prepare cards

Dry blood spot (DBS) cards arrive and are checked for completeness.



Batching & scanning

Different card types are organized for processing and scanned.




Data transfer

Revvity Instrument Hub transfers data to Transcribe AI in the cloud.



Digital conversion

The handwritten sections are converted into a digital format via Transcribe AI.



Data transfer

The Revvity Instrument Hub seamlessly transfers data between Revvity Transcribe AI and your lab system.



Validate results

A quick human review ensures accuracy before the data is finalized.

Automatic

Elevate your lab's productivity with Revvity Transcribe AI

Why use Revvity Transcribe AI

- For use with handwritten text on Test request forms (TRF) including DBS cards.
- ISO27001 compliant for secure cloud-based operations.
- Reduce the need for dedicated staff by automating the transcription.

Security and privacy

- Data is not stored in cloud. The images are read utilizing cloud service but images are not retained in cloud or used in training of models.
- Revvity Instrument Hub, Revvity Transcribe AI application, and the cloud environment are ISO27001 compliant.

Order Revvity Transcribe AI

| Part number | Description |
|-------------|------------------------|
| 5024-1000 | Revvity Instrument Hub |
| 5024-2000 | Transcribe AI |

Technical requirements

Scanner:

- **Image size:** 1500px x 1000px or larger
- **Resolution:** 96 DPI or higher
- **File size:** Under 5MB (if larger, the system compresses them)
- **File formats:** TIFF, JPEG, or JPG

Revvity Instrument Hub is needed to use Revvity Transcribe AI

