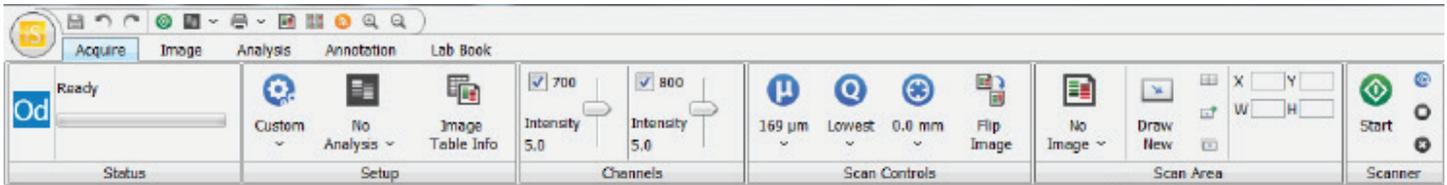




# Image Studio 4.0 Imaging Software

## Image Acquisition and Import

- **Application-Driven Ribbon Interface (Figure 1)**
  - Each analysis type has a corresponding ribbon with tools for analysis and formatting that are only relevant for that application



**Figure 1.** Image Studio is a ribbon-based application that displays analysis and formatting tools for user selection and implementation.

- **Ability to scan multiple areas in a single scan**
  - With an enhanced acquisition interface, the user can divide a single scan into multiple scan areas
- **Ability to import .tif, .jpg and .png files**

## Data Analysis

- **Definition of Lanes and Bands**
  - Simply define the sample area and the total number of lanes
- **Placement of Shapes by Simple Point-and-Click Method**
  - Click on a band to see a rectangle placed around it automatically
- **Rotation of Objects**
  - If a band or blot are not in the desired orientation, objects can be rotated as a group or individually
- **Data Analysis Tools**
  - The user has full control of the data display and can rearrange each column through sorting and filtering
  - Create customized charts
- **Annotation Controls**
  - Simple to change fonts and controls, including rotational control of text and arrows
- **Linking of Look-Up-Tables**
  - View different scans under identical display conditions
- **Supports many types of analysis, including:**
  - Western Blots
  - In-Cell Western™ Assays
  - Multi-Well Plates
  - Coomassie Gel Documentation
  - MPX™ Western Blots
  - EMSAs
  - Tissue Imaging
  - Grid Arrays
  - DNA Gels
  - Small Animal Imaging

