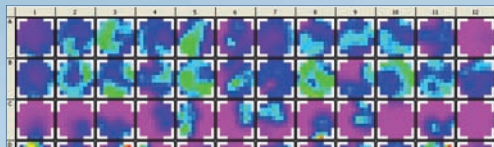


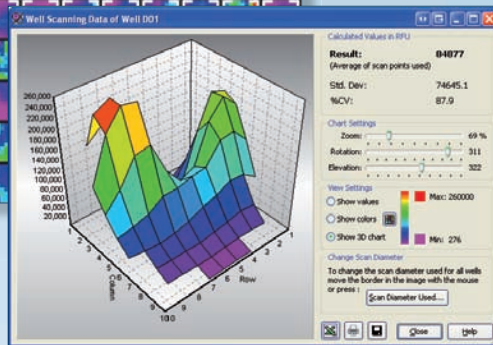
## PERFORM NEW ASSAYS FASTER AND EASIER DATA ANALYSIS SOFTWARE

### 3D Well Mapping

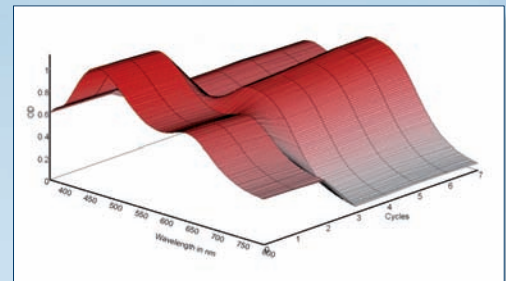


Well scanning of non-confluent cells

3D well map of non-confluent cells in a 96-well plate



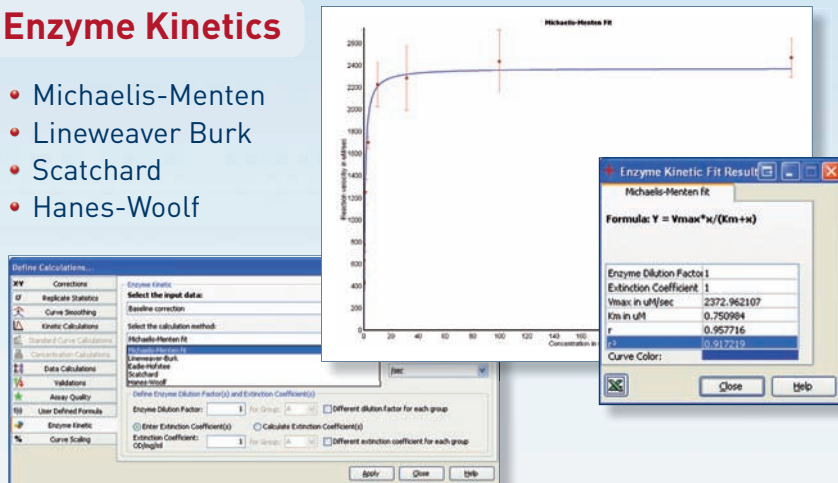
Full-spectrum kinetic absorbance data for an enzymatic assay



3D well map of a Bradford absorbance spectrum taken over 60 seconds – 475nm peak decreases and 595nm peak increases

### Enzyme Kinetics

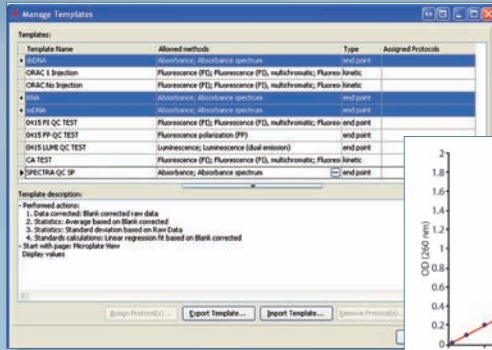
- Michaelis-Menten
- Lineweaver Burk
- Scatchard
- Hanes-Woolf



Michaelis-Menten calculation produces a graph,  $K_m$ , and  $V_{max}$  values for enzymatic assays

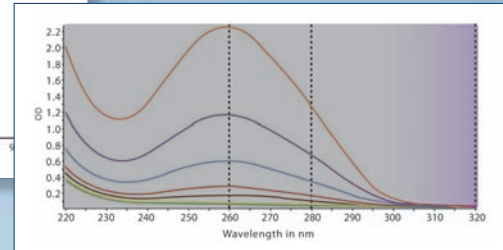
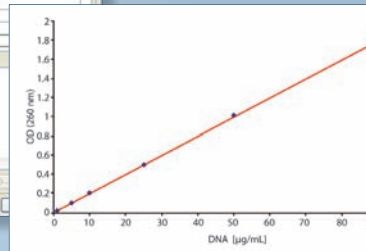


## Assay Calculation Templates



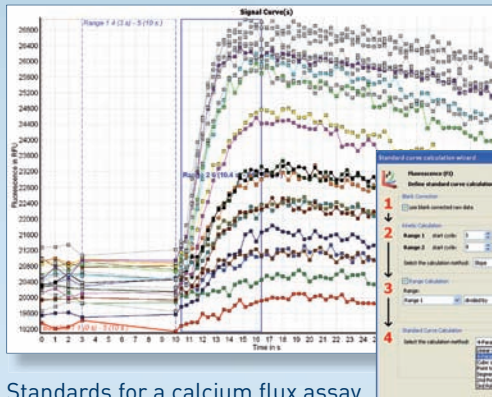
Automatic DNA/RNA concentration calculation

Determines DNA concentration from absorbance 260/280 readings (<0.1ug/mL at 3 uL in 384-well plates)

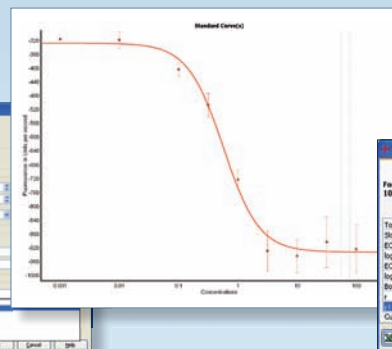


DNA calculation template uses a preset standard curve or creates a new one from user's data to determine DNA concentration of unknowns.

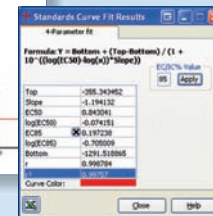
## Standard Curve Wizard



Calculates a curve from known standards using a wizard

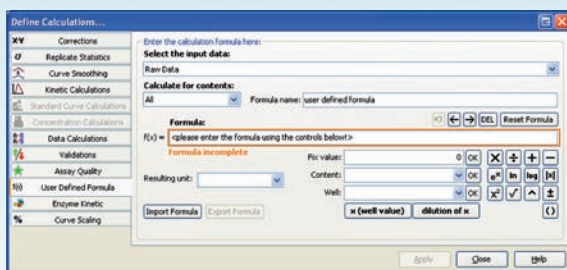


Wizard helps create a 4-parameter fit;  $r^2$ ,  $EC_{50}$  and  $EC_{85}$  values are determined

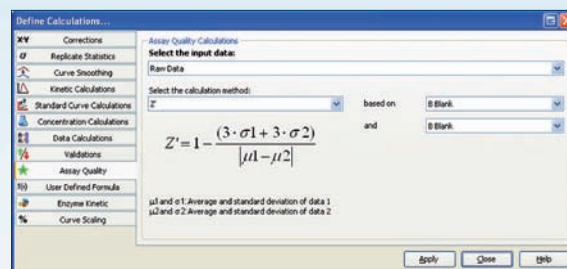


Standards for a calcium flux assay

## User-Defined and Standard Formulas for Calculations



Z' is a standard calculation in MARS that is used to evaluate HTS assays.



The MARS formula generator allows users to define and perform any desired calculation

## Additional Features

- Full Spectrum Absorbance Measurements
- Signal-to-noise, signal-to-blank, and F% calculations
- Determine any EC/IC value
- Simple export to Excel and ASCII
- Print to PDFs
- Easy layout edit
- 21 CFR part 11 compliant