

## In Silico Prediction

# chemTarget™ - Predicts Biological Target Interaction Directly from Chemical Structure

## Background Information



- Predicts binding affinity, inhibition constants or other measures of interaction with biological targets, directly from chemical structure.
- Uses Cyprotex's unique pattern recognition software to build models from existing data sets (provided by the customer or from the literature).
- Analyses approximately 10,000 descriptors using linear, random forest, neural network and nearest neighbour methods.
- Provides clinically relevant binding/inhibition/activation when used in combination with the pharmacokinetic predictor, chemPK™.
- Provides an early-stage filter for directing chemistry and prioritising screening.

### chemTarget™ Input Requirements

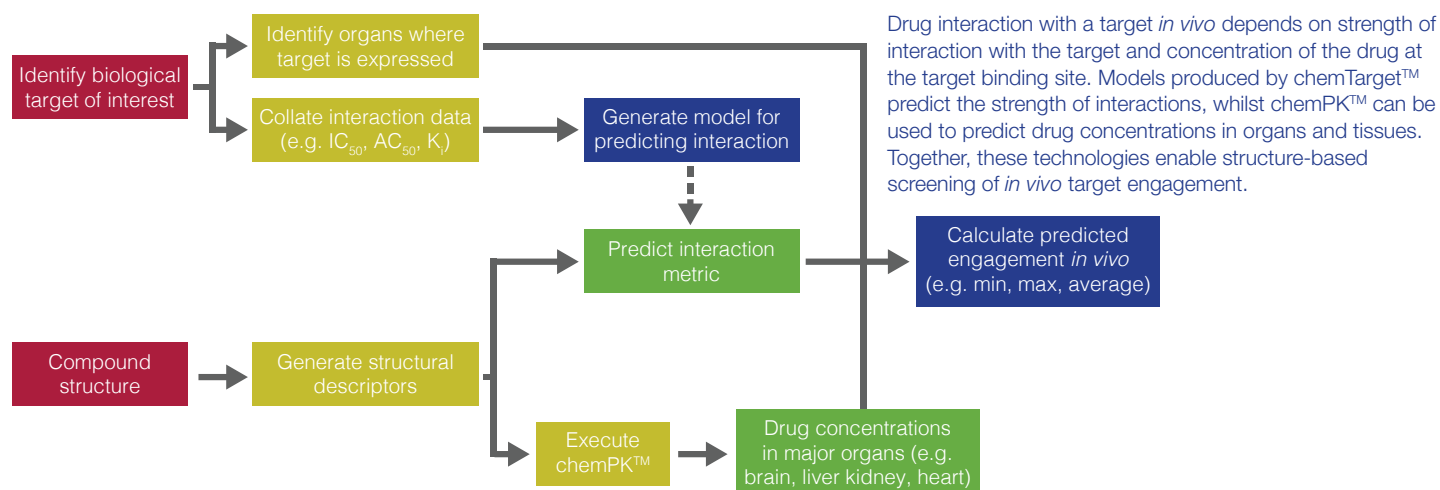
- Chemical structure, e.g., SMILES, mol or sdf.
- Location(s) of *in vivo* target expression.
- Existing target interaction data (e.g.  $IC_{50}$ ,  $AC_{50}$  or  $K_i$ ).

### chemTarget™ Data Delivery

- Predicted target interaction.
- Predicted engagement *in vivo* for specified dose-regimen(s) (minimum, maximum, average) if used in combination with chemPK™.

**Figure 1**

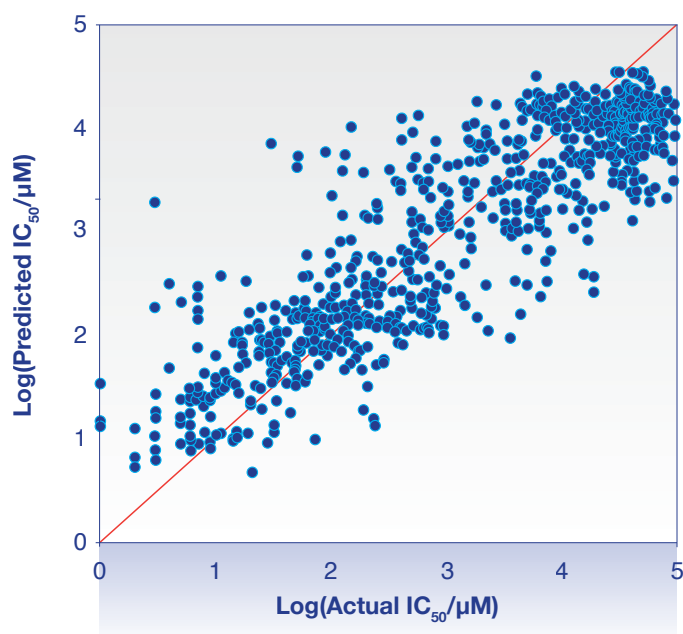
Schematic illustrating how chemTarget™ can be integrated with chemPK™ to predict clinically relevant biological target interaction.



## Performance of chemTarget™ predictions

**Figure 2**

Prediction of JNK3 binding affinity from chemical structure. Results are from 10 repeats of 10-fold cross-validation for a set of 697 compounds.



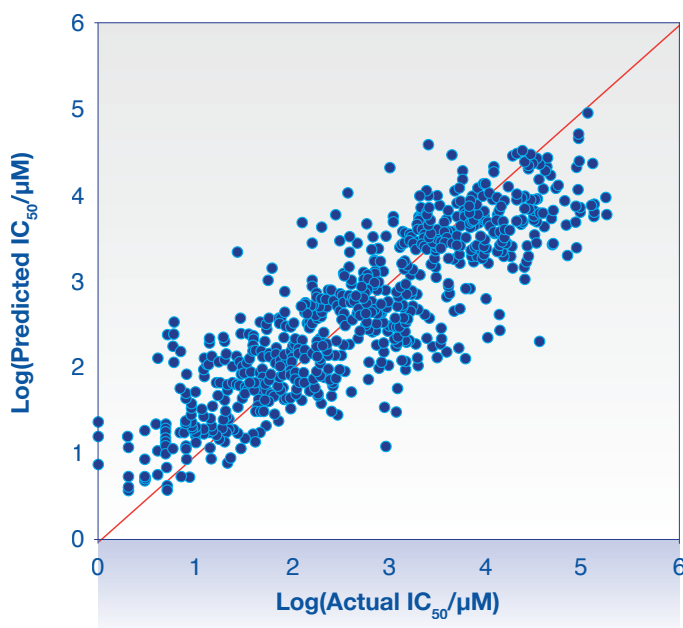
RMSE*	0.68
R <sup>2</sup>	0.76
Spearman rank correlation coefficient	0.85

\*RMSE = root mean square error

JNK3 is a potential therapeutic target for several neurodegenerative disorders. chemTarget™ predicts JNK3 inhibition directly from structure with a repeated cross-validation R<sup>2</sup> of 0.76 for a set of 697 compounds

**Figure 3**

Prediction of MK2 binding affinity from chemical structure. Results are from 10 repeats of 10-fold cross-validation for a set of 670 compounds.



RMSE*	0.64
R <sup>2</sup>	0.72
Spearman rank correlation coefficient	0.85

\*RMSE = root mean square error

MK2 (mitogen-activated protein kinase (MAPK)-activated protein kinase 2) is a potential therapeutic target in inflammatory disease. chemTarget™ predicts MK2 inhibition directly from structure with a repeated cross-validation R<sup>2</sup> of 0.72 for a set of 670 compounds.

Contact [enquiries@cyprotex.com](mailto:enquiries@cyprotex.com) to discuss your project.