

“Rational” solutions for 4 key issues in protein crystallography

