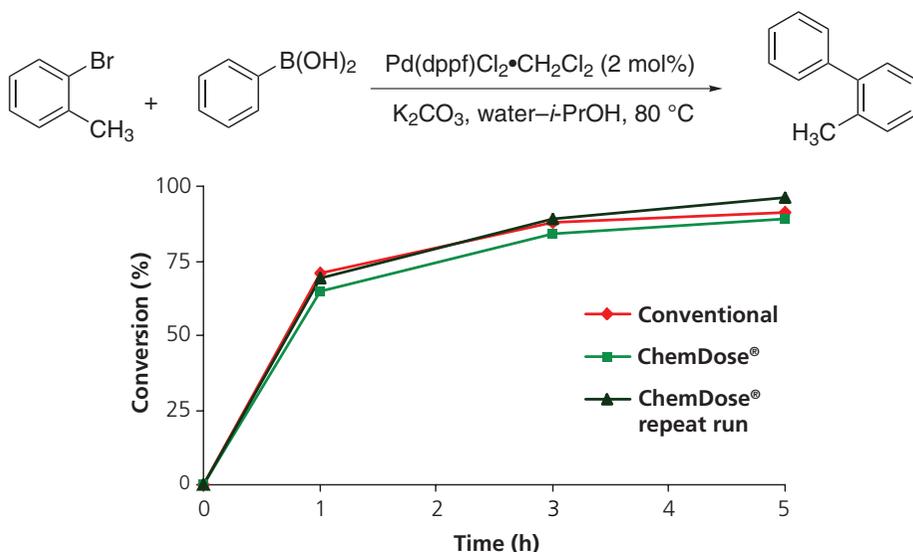


Frequently Asked Questions (FAQs)

How do ChemDose® tablets compare to conventional catalysts, ligands or reagents?

As illustrated in the graph below (a Suzuki coupling reaction), typically the kinetics and conversions are quite similar.



What are the ChemDose® tablets made from?

How big are the tablets?

The tablets themselves are composed of a chemically inert magnesium aluminosilicate matrix, along with the adsorbed reagent or catalyst. All tablets are approximately 5 millimeters in diameter.

Where do I find information on the exact loading for a given lot of a ChemDose® product?

The certificate of analysis (C of A) contains lot specific loading information, along with the method used for the loading determination.

How should I dispose of spent ChemDose® tablets?

They should be treated as solid waste, contaminated with reaction species. The matrix is non-toxic and environmentally benign.

How should I store ChemDose® tablets?

As with most catalysts and ligands, as a general precaution it is best to store them in a refrigerator and/or under an inert atmosphere to maintain activity. Some catalysts, such as Pd(dppf)Cl₂·CH₂Cl₂, are more robust and can be handled in air without issue. Each ChemDose® label provides specific handling information.

Can I use them in any solvent?

Yes, including aqueous solvents.

Do the tablets disintegrate during the course of the reaction?

They are normally recovered intact, however, they may break apart somewhat in aqueous reactions. Any disintegrated tablets can be removed from the reaction via filtration.

What is the pH stability of the ChemDose® tablets?

The ChemDose® tablets are mildly basic and will dissolve in the presence of mineral acids.

Can I use them in a microwave or pressure vessel?

Absolutely. Please see sigma-aldrich.com/chemdose for examples of cross-coupling reactions involving the use of microwaves.



SIGMA-ALDRICH