

**Application Guide for Palladium
Catalyzed Cross-Coupling Reactions**

Add Aldrich



Common Cross-Coupling Reactions

Negishi (C-C)



R¹ = alkenyl, (hetero)aryl, alkynyl, acyl
 R² = alkenyl, (hetero)aryl, alkynyl, alkyl
 X = Cl, Br, I, OTf, OAc

Heck (C-C)



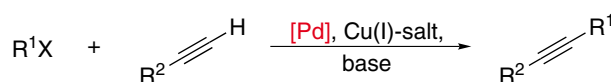
R¹ = alkenyl, (hetero)aryl, alkyl (w/o β-hydrogens)
 R² = alkenyl, (hetero)aryl, alkyl
 X = Cl, Br, I, OTf, OTs, N₂⁺

Suzuki (C-C)



R¹ = alkenyl, (hetero)aryl, alkyl
 R² = alkenyl, (hetero)aryl, alkynyl, alkyl
 Y = alkyl, OH, O-alkyl
 X = Cl, Br, I, OTf

Sonogashira (C-C)



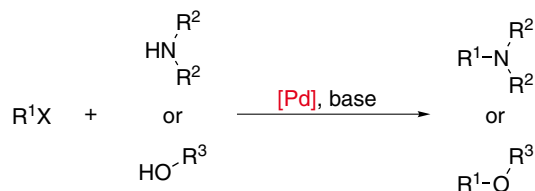
R¹ = alkenyl, (hetero)aryl
 R² = H, alkenyl, (hetero)aryl, alkyl, SiR₃
 X = Cl, Br, I, OTf

Stille (C-C)



R¹ = alkenyl, (hetero)aryl, acyl
 R² = alkenyl, (hetero)aryl, alkyl
 X = Cl, Br, I, OTf

Buchwald-Hartwig (C-N and/or C-O)


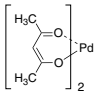

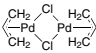

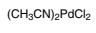

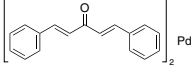

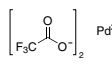

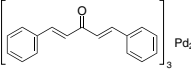
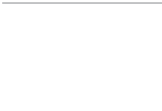
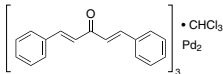

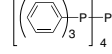

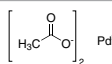

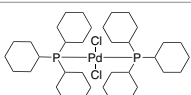

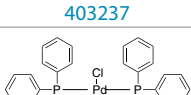


R¹ = (hetero)aryl
 R² = 1° or 2° aryl or alkyl
 R³ = 1°, 2°, or 3° alkyl, aryl
 X = Cl, Br, I, OTf


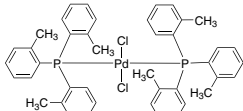









References:

- Negishi, E., *Handbook of Organopalladium Chemistry for Organic Synthesis*; John Wiley & Sons: New York, NY, 2002.
- Tsuji, J., *Palladium Reagents and Catalysts: New Perspectives for the 21st Century*; John Wiley & Sons: Chichester, West Sussex, England, 2004.
- Kürti, L. and Czakó, B. *Strategic Applications of Named Reactions in Organic Synthesis*; Elsevier Academic Press: Burlington, MA, 2005.
- Li, J. J. *Named Reactions for Homologations*; John Wiley & Sons: Hoboken, NJ, 2009.


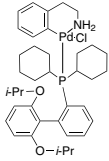

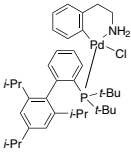

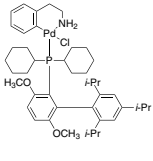

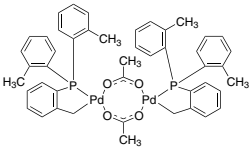

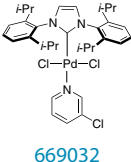
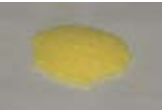
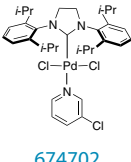

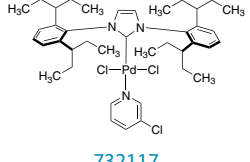
Palladium Catalysts and Precursors

Structure and Cat. No.	Common Name	CAS No.	Mol. Wt.	Color	Stability	Negishi	Suzuki	Stille	Heck	Sonogashira	Buchwald-Hartwig
  209015	Pd(acac) ₂	14024-61-4	304.64	Yellow	Air Stable		X		X		X
  222380	[Pd(allyl)Cl] ₂	12012-95-2	365.89	Yellow	Air/Moisture Sensitive		X	X	X	X	
  225657	Pd(MeCN) ₂ Cl ₂	14592-56-4	259.43	Dark Orange	Air Stable	X	X	X		X	
  227994	Pd(dba) ₂	32005-36-0	575.00	Black	Air/Moisture Sensitive	X	X	X	X		X
  299685	Pd(TFA) ₂	42196-31-6	332.45	Brown	Air Stable		X		X	X	X
  328774	Pd ₂ (dba) ₃	51364-51-3	915.72	Dark Purple to Black	Air Stable	X	X		X		X
  366315 (chloroform adduct of 328774)	Pd ₂ (dba) ₃ · CHCl ₃	52522-40-4	1035.10	Purple	Air Stable		X		X		X
  216666 697265	Pd(PPh ₃) ₄	14221-01-3	1155.56	Yellow	Air/Moisture Sensitive	X	X	X		X	X
  520764	Pd(OAc) ₂	3375-31-3	224.51	Orange-Brown	Air Stable	X	X	X	X		X
  403237	Pd(PCy ₃) ₂ Cl ₂	29934-17-6	738.18	Yellow	Air Stable		X		X	X	X
  412740	Pd(PPh ₃) ₂ Cl ₂	13965-03-2	701.90	Yellow	Air/Moisture Sensitive	X	X	X	X	X	

Application Guide for Palladium Catalyzed Cross-Coupling Reactions

Structure and Cat. No.	Common Name	CAS No.	Mol. Wt.	Color	Stability	Negishi	Suzuki	Stille	Heck	Sonogashira	Buchwald-Hartwig
  453048	Pd[P(o-tol) ₃] ₂ Cl ₂	40691-33-6	786.06	Yellow	Air/Moisture Sensitive	X	X	X	X		X
 678740	Pd(amphos)Cl ₂	887919-35-9	708.07	Yellow	Air Stable	X	X				
 697230	Pd(dppf)Cl ₂	72287-26-4	731.70	Red	Air Stable	X	X	X			
 379670	Pd(dppf)Cl ₂ · CH ₂ Cl ₂	95464-05-4	816.64	Red	Air Stable	X	X	X			
 701602	Pd(dtbpf)Cl ₂	95408-45-0	651.75	Brown	Air Stable		X		X		
 339768	Pd(MeCN) ₄ (BF ₄) ₂	21797-13-7	444.24	Dark Yellow	Air/Moisture Sensitive	X	X		X		X
 205877	PdBr ₂	13444-94-5	266.23	Dark Brown	Air Stable				X	X	X
 520659	PdCl ₂	7647-10-1	177.33	Red-Brown	Air Stable		X	X	X	X	
 NEW 704946	(SPhos) palladium(II) phenethylamine chloride (1:1 MTBE solvate)	N/A	760.72	White	Air-Stable						X
 NEW 704954	(XPhos) palladium(II) phenethylamine chloride	1028206-56-5	738.76	White	Air-Stable						X

Palladium Catalysts and Precursors

	Structure and Cat. No.	Common Name	CAS No.	Mol. Wt.	Color	Stability	Negishi	Suzuki	Stille	Heck	Sonogashira	Buchwald-Hartwig
 NEW	 707589	(RuPhos) palladium(II) phenethylamine chloride	1028206-60-1	728.68	White	Air-Stable						X
 NEW	 708739	(t-BuXPhos) palladium(II) phenethylamine chloride	1142811-12-8	686.69	White	Air-Stable						X
 NEW	 718750	(BrettPhos) palladium(II) phenethylamine chloride	1148148-01-9	798.81	White	Air Stable						X
	 671916	Herrmann's Catalyst or cataCXium® C	172418-32-5	937.64	Light yellow	Air Stable				X		
	 669032	Pd-PEPPSI™-IPr	905459-27-0	679.46	Yellow	Air Stable	X	X			X	X
	 674702	Pd-PEPPSI™-SIPr	927706-57-8	681.47	Yellow	Air Stable	X	X				X
 NEW	 732117	Pd-PEPPSI™-IPent	1158652-41-5	791.67	Yellow	Air Stable	X	X	X			

Sigma-Aldrich® Worldwide Offices

Argentina

Free Tel: 0810 888 7446
Tel: (+54) 11 4556 1472
Fax: (+54) 11 4552 1698

Australia

Free Tel: 1800 800 097
Free Fax: 1800 800 096
Tel: (+61) 2 9841 0555
Fax: (+61) 2 9841 0500

Austria

Tel: (+43) 1 605 81 10
Fax: (+43) 1 605 81 20

Belgium

Free Tel: 0800 14747
Free Fax: 0800 14745
Tel: (+32) 3 899 13 01
Fax: (+32) 3 899 13 11

Brazil

Free Tel: 0800 701 7425
Tel: (+55) 11 3732 3100
Fax: (+55) 11 5522 9895

Canada

Free Tel: 1800 565 1400
Free Fax: 1800 265 3858
Tel: (+1) 905 829 9500
Fax: (+1) 905 829 9292

Chile

Tel: (+56) 2 495 7395
Fax: (+56) 2 495 7396

China

Free Tel: 800 819 3336
Tel: (+86) 21 6141 5566
Fax: (+86) 21 6141 5567

Czech Republic

Tel: (+420) 246 003 200
Fax: (+420) 246 003 291

Denmark

Tel: (+45) 43 56 59 00
Fax: (+45) 43 56 59 05

Finland

Tel: (+358) 9 350 9250
Fax: (+358) 9 350 92555

France

Free Tel: 0800 211 408
Free Fax: 0800 031 052
Tel: (+33) 474 82 28 88
Fax: (+33) 474 95 68 08

Germany

Free Tel: 0800 51 55 000
Free Fax: 0800 64 90 000
Tel: (+49) 89 6513 0
Fax: (+49) 89 6513 1160

Hungary

Ingyenes telefonszám: 06 80 355 355
Ingyenes fax szám: 06 80 344 344
Tel: (+36) 1 235 9063
Fax: (+36) 1 269 6470

India

Telephone

Bangalore: (+91) 80 6621 9400
New Delhi: (+91) 11 4358 8000
Mumbai: (+91) 22 2570 2364
Hyderabad: (+91) 40 4015 5488
Kolkata: (+91) 33 4013 8003

Fax

Bangalore: (+91) 80 6621 9550
New Delhi: (+91) 11 4358 8001
Mumbai: (+91) 22 4087 2364
Hyderabad: (+91) 40 4015 5488
Kolkata: (+91) 33 4013 8000

Ireland

Free Tel: 1800 200 888
Free Fax: 1800 600 222
Tel: (+353) 402 20370
Fax: (+353) 402 20375

Israel

Free Tel: 1 800 70 2222
Tel: (+972) 8 948 4100
Fax: (+972) 8 948 4200

Italy

Free Tel: 800 827 018
Tel: (+39) 02 3341 7310
Fax: (+39) 02 3801 0737

Japan

Tel: (+81) 3 5796 7300
Fax: (+81) 3 5796 7315

Korea

Free Tel: (+82) 80 023 7111
Free Fax: (+82) 80 023 8111
Tel: (+82) 31 329 9000
Fax: (+82) 31 329 9090

Malaysia

Tel: (+60) 3 5635 3321
Fax: (+60) 3 5635 4116

Mexico

Free Tel: 01 800 007 5300
Free Fax: 01 800 712 9920
Tel: (+52) 722 276 1600
Fax: (+52) 722 276 1601

The Netherlands

Free Tel: 0800 022 9088
Free Fax: 0800 022 9089
Tel: (+31) 78 620 5411
Fax: (+31) 78 620 5421

New Zealand

Free Tel: 0800 936 666
Free Fax: 0800 937 777
Tel: (+61) 2 9841 0555
Fax: (+61) 2 9841 0500

Norway

Tel: (+47) 23 17 60 00
Fax: (+47) 23 17 60 10

Poland

Tel: (+48) 61 829 01 00
Fax: (+48) 61 829 01 20

Portugal

Free Tel: 800 202 180
Free Fax: 800 202 178
Tel: (+351) 21 924 2555
Fax: (+351) 21 924 2610

Russia

Tel: (+7) 495 621 5828
Fax: (+7) 495 621 6037

Singapore

Tel: (+65) 6779 1200
Fax: (+65) 6779 1822

Slovakia

Tel: (+421) 255 571 562
Fax: (+421) 255 571 564

South Africa

Free Tel: 0800 1100 75
Free Fax: 0800 1100 79
Tel: (+27) 11 979 1188
Fax: (+27) 11 979 1119

Spain

Free Tel: 900 101 376
Free Fax: 900 102 028
Tel: (+34) 91 661 99 77
Fax: (+34) 91 661 96 42

Sweden

Tel: (+46) 8 742 4200
Fax: (+46) 8 742 4243

Switzerland

Free Tel: 0800 80 00 80
Free Fax: 0800 80 00 81
Tel: (+41) 81 755 2828
Fax: (+41) 81 755 2815

United Kingdom

Free Tel: 0800 717 181
Free Fax: 0800 378 785
Tel: (+44) 1747 833 000
Fax: (+44) 1747 833 313

United States

Toll-Free: 800 325 3010
Toll-Free Fax: 800 325 5052
Tel: (+1) 314 771 5765
Fax: (+1) 314 771 5757

Vietnam

Tel: (+84) 3516 2810
Fax: (+84) 6258 4238

Internet

sigma-aldrich.com

*Enabling Science to
Improve the Quality of Life*

Order/Customer Service (800) 325-3010 • Fax (800) 325-5052
Technical Service (800) 325-5832 • sigma-aldrich.com/techservice
Development/Custom Manufacturing Inquiries **SAFC® (800) 244-1173**
Safety-related Information sigma-aldrich.com/safetycenter

World Headquarters
3050 Spruce St.
St. Louis, MO 63103
(314) 771-5765
sigma-aldrich.com

©2011 Sigma-Aldrich Co. All rights reserved. SIGMA, SAFC, SIGMA-ALDRICH, ALDRICH and SUPELCO are registered trademarks of Sigma-Aldrich Biotechnology LP, an affiliate of Sigma-Aldrich Co., FLUKA is a registered trademark of Fluka GmbH. Sigma brand products are sold through Sigma-Aldrich, Inc. Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip. cataCXium is a registered trademark of Evonik Degussa GmbH. PEPPSI is a trademark of Total Synthesis Ltd.

NKO
75683-510273
1021