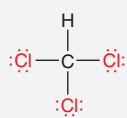
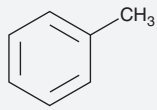
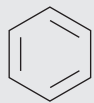
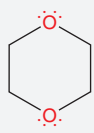
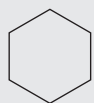
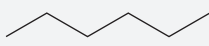
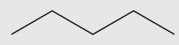


COMMON NONPOLAR SOLVENTS

Nonpolar solvents have low dielectric constants and are poor solvents for charged species. Diethyl ether, used for Grignard reactions, is one notable exception since its lone pairs can help to solvate the Mg cation.

Solvent	Structure	Dielectric Constant	Dipole Moment	Boiling Point (°C)
Chloroform		4.81	1.15	61.7
Diethyl ether	$\text{H}_3\text{CH}_2\text{C}-\ddot{\text{O}}-\text{CH}_2\text{CH}_3$	4.33	1.15	34.6
Toluene		2.38	0.375	110.6
Benzene		2.28	0	80.1
1,4-Dioxane		2.25	0.45	102
Cyclohexane		2.02	0	80.7
Hexane		1.88	1.08	69
Pentane		1.84	0.007	36.1