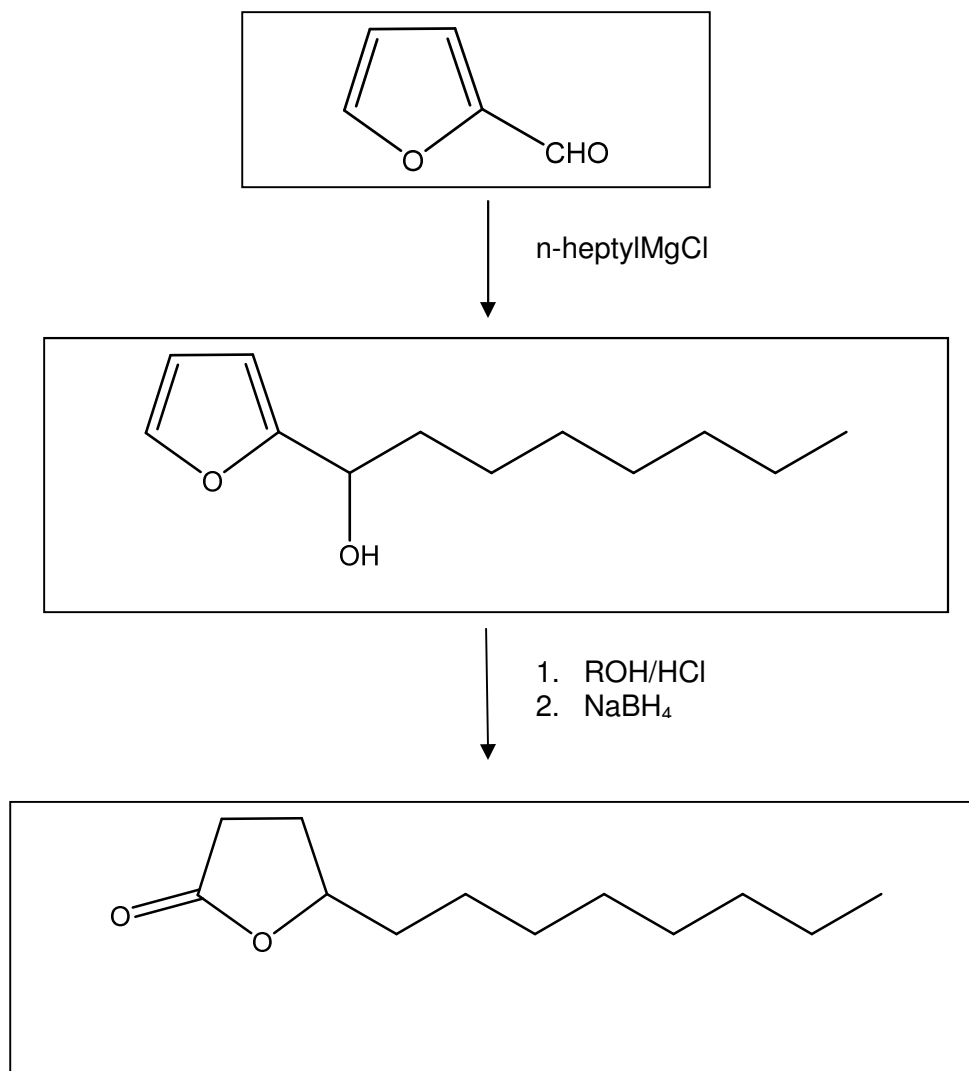


Process Flow

The process flow for the following product is listed below.

W240001 γ -Dodecalactone, $\geq 97\%$, FCC, FG



Sigma-Aldrich warrants that, as of the above date, this product conformed to the information contained in this publication. For further inquiries, please contact Technical Service. Purchaser must determine the suitability of the product for its particular use. See sigma-aldrich.com, the reverse side of invoice or packing slip for additional terms and conditions of sale.